

# Accessibility Conformance Report

## Revised Section 508 Edition

### (Based on VPAT® Version 2.5)

**Name of Product/Version:** Belysa® analysis software version 1.2.2

**Report Date:** 4/3/2024

**Product Description:** Belysa® analysis software displays single and multiplex data that allows the user to immediately visualize their standard curve and the distribution of standard points, controls, and samples.

**Contact Information:** [ImmunoassayTS@milliporesigma.com](mailto:ImmunoassayTS@milliporesigma.com)

**Notes:** Evaluated on Windows 10 platform with Narrator

**Evaluation Methods Used:** Testing is based on knowledge of general product functionality and general knowledge of accessibility technology.

#### Applicable Standards/Guidelines

This report covers the degree of conformance for the following accessibility standard/guidelines:

Standard/Guideline	Included In Report
<a href="#">Web Content Accessibility Guidelines 2.0</a>	Level A (Yes) Level AA (Yes) Level AAA (No)
<a href="#">Revised Section 508 standards published January 18, 2017 and corrected January 22, 2018</a>	(Yes)

## Terms

The terms used in the Conformance Level information are defined as follows:

- **Supports:** The functionality of the product has at least one method that meets the criterion without known defects or meets with equivalent facilitation.
- **Partially Supports:** Some functionality of the product does not meet the criterion.
- **Does Not Support:** The majority of product functionality does not meet the criterion.
- **Not Applicable:** The criterion is not relevant to the product.
- **Not Evaluated:** The product has not been evaluated against the criterion. This can only be used in WCAG Level AAA criteria.

## WCAG 2.0 Report

Tables 1 and 2 also document conformance with Revised Section 508:

- Chapter 5 – 501.1 Scope, 504.2 Content Creation or Editing
- Chapter 6 – 602.3 Electronic Support Documentation

Note: When reporting on conformance with the WCAG 2.0 Success Criteria, they are scoped for full pages, complete processes, and accessibility-supported ways of using technology as documented in the [WCAG 2.0 Conformance Requirements](#).

**Table 1: Success Criteria, Level A**

Notes: Belysa software application itself is considered "Software" for the purposes of Conformance Level labeling. Integrated browser-based user documentation has been classified as "Electronic Docs". A User Guide is also available in PDF format.

Criteria	Conformance Level	Remarks and Explanations
<b><u>1.1.1 Non-text Content</u></b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"><li>• 501 (Web)(Software)</li><li>• 504.2 (Authoring Tool)</li><li>• 602.3 (Support Docs)</li></ul>	Electronic Docs: <b>Partially Supports</b> Software: <b>Supports</b>	Electronic Docs: Many images in documentation are presented next to text explanations, but some screenshots are not captioned/do not have alternative text  Software: Graphical data presented in equivalent tabular form; controls are labeled with text that describes its purpose
<b><u>1.2.1 Audio-only and Video-only (Prerecorded)</u></b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"><li>• 501 (Web)(Software)</li><li>• 504.2 (Authoring Tool)</li><li>• 602.3 (Support Docs)</li></ul>	Electronic Docs: <b>Not Applicable</b> Software: <b>Not Applicable</b>	Electronic Docs: No audio-only or video-only content  Software: No audio-only or video-only content
<b><u>1.2.2 Captions (Prerecorded)</u></b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"><li>• 501 (Web)(Software)</li><li>• 504.2 (Authoring Tool)</li><li>• 602.3 (Support Docs)</li></ul>	Electronic Docs: <b>Not Applicable</b> Software: <b>Not Applicable</b>	Electronic Docs: No prerecorded audio or synchronized media  Software: No prerecorded audio or synchronized media Authoring Tool:
<b><u>1.2.3 Audio Description or Media Alternative (Prerecorded)</u></b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"><li>• 501 (Web)(Software)</li><li>• 504.2 (Authoring Tool)</li><li>• 602.3 (Support Docs)</li></ul>	Electronic Docs: <b>Not Applicable</b> Software: <b>Not Applicable</b>	Electronic Docs: No time-based media or prerecorded video  Software: No time-based media or prerecorded video
<b><u>1.3.1 Info and Relationships</u></b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"><li>• 501 (Web)(Software)</li><li>• 504.2 (Authoring Tool)</li><li>• 602.3 (Support Docs)</li></ul>	Electronic Docs: <b>Partially Supports</b> Software: <b>Partially Supports</b>	Electronic Docs: Content rendered using semantic HTML, though not all accessibility features supported.  Software: Application is a compiled Windows native

Criteria	Conformance Level	Remarks and Explanations
		program. Windows 10 Navigator able to determine some menu labels, menu state, table rows/columns, and other aspects of information displayed on screen, but not all functions compatible.
<u><b>1.3.2 Meaningful Sequence</b></u> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b> Software: <b>Supports</b>	Electronic Docs: CSS layouts used, tables not used for positioning  Software: In general, the software does not employ content whose meaning is significantly dependent on specific sequence
<u><b>1.3.3 Sensory Characteristics</b></u> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b> Software: <b>Supports</b>	Electronic Docs: Does not contain content that relies solely upon sensory characteristics of components  Software: Graphical content such as well group designations (Standard, Control, Unknown) and well message labels ("EXT", "ND") are presented using symbols that differ in both shape and color, and with accompanying text
<u><b>1.4.1 Use of Color</b></u> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b> Software: <b>Supports</b>	Electronic Docs: Color not used as the only means of conveying information.  Software: Color is used to differentiate types of wells (Standard, Control, Unknown), but shape and text labeling are also differentiated in this case
<u><b>1.4.2 Audio Control</b></u> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Not Applicable</b> Software: <b>Not Applicable</b>	Electronic Docs: No audio content Software: No audio content

Criteria	Conformance Level	Remarks and Explanations
<b>2.1.1 Keyboard</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b> Software: <b>Partially Supports</b>	Electronic Docs: All functionality accessible via keyboard  Software: Keyboard shortcuts are provided for many common operations (opening/closing files, changing tabs, editing information), but other operations (adding groups to experiments) are missing keyboard functionality
<b>2.1.2 No Keyboard Trap</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b> Software: <b>Does Not Support</b>	Electronic Docs: No keyboard traps, standard web browser interface  Software: Keyboard traps were observed in the software and required non-standard arrow/tab keys to escape
<b>2.2.1 Timing Adjustable</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b> Software: <b>Supports</b>	Electronic Docs: No functions with time limits Software: No functions with time limits
<b>2.2.2 Pause, Stop, Hide</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b> Software: <b>Supports</b>	Electronic Docs: No moving, blinking, or auto-updating content Software: No moving, blinking, or auto-updating content
<b>2.3.1 Three Flashes or Below Threshold</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b> Software: <b>Supports</b>	Electronic Docs: No flashing content Software: No flashing content

Criteria	Conformance Level	Remarks and Explanations
<b>2.4.1 Bypass Blocks</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>501 (Web)(Software) – Does not apply to non-web software</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs) – Does not apply to non-web docs</li> </ul>	Electronic Docs: <b>Supports</b>	Electronic Docs:
<b>2.4.2 Page Titled</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>501 (Web)(Software)</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b> Software: <b>Supports</b>	Electronic Docs: Pages display title in title element Software: All page tabs are labeled with section title ("Results", "Raw Data", etc.)
<b>2.4.3 Focus Order</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>501 (Web)(Software)</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b> Software: <b>Does Not Support</b>	Electronic Docs: Navigation sequence does not affect meaning or operation  Software: Focus order does not preserve optimal operability in many areas
<b>2.4.4 Link Purpose (In Context)</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>501 (Web)(Software)</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b> Software: <b>Supports</b>	Electronic Docs: Links generally reflect topic or section to which they will navigate when clicked  Software: No use of text links other than About screen, where purpose is labeled
<b>3.1.1 Language of Page</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>501 (Web)(Software)</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b> Software: <b>Does Not Support</b>	Electronic Docs: HTML labeled as "en-US" Software: Application is not HTML based and does not provide an explicit API for determining language
<b>3.2.1 On Focus</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>501 (Web)(Software)</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b>  Software: <b>Supports</b>	Electronic Docs: No change of context on focus  Software: No change of context on focus

Criteria	Conformance Level	Remarks and Explanations
<b>3.2.2 On Input</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b>  Software: <b>Supports</b>	Electronic Docs: No change of context on input  Software: No change of context on input
<b>3.3.1 Error Identification</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b>  Software: <b>Partially Supports</b>	Electronic Docs: No input required.  Software: In many cases invalid input is prevented from being entered (e.g., prevent letters in numeric inputs). When invalid input is able to be entered and detected, typically the control reverts to its last valid value and does not display an error.
<b>3.3.2 Labels or Instructions</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b>  Software: <b>Does Not Support</b>	Electronic Docs: No user input required.  Software: Labels and instructions typically not provided for user input. User must rely on documentation or context.
<b>4.1.1 Parsing</b> (Level A) Also applies to: WCAG 2.0 – Always answer 'Supports' Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	<b>Supports</b>	For WCAG 2.0 and the 508 standards, the September 2023 errata update indicates this criterion is always supported. See the <a href="#">WCAG 2.0 Editorial Errata</a> .
<b>4.1.2 Name, Role, Value</b> (Level A) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b>  Software: <b>Does Not Support</b>	Electronic Docs: Minimal/no use of custom user interface components  Software: User interface components generally not able to provide state or role information to assistive technologies



**Table 2: Success Criteria, Level AA**

Notes:

Criteria	Conformance Level	Remarks and Explanations
<b><u>1.2.4 Captions (Live)</u></b> (Level AA) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>501 (Web)(Software)</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Not Applicable</b>  Software: <b>Not Applicable</b>	Electronic Docs: No live audio content  Software: No live audio content
<b><u>1.2.5 Audio Description (Prerecorded)</u></b> (Level AA) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>501 (Web)(Software)</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Not Applicable</b>  Software: <b>Not Applicable</b>	Electronic Docs: No prerecorded video content  Software: No prerecorded video content
<b><u>1.4.3 Contrast (Minimum)</u></b> (Level AA) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>501 (Web)(Software)</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Partially Supports</b>  Software: <b>Partially Supports</b>	Electronic Docs: Text is generally black on white background, but no specific contrast controls exist Software: Text is generally black on white background, dark mode option exists, but no specific contrast controls exist
<b><u>1.4.4 Resize text</u></b> (Level AA) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>501 (Web)(Software)</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Does Not Support</b>  Software: <b>Does Not Support</b>	Electronic Docs: No independent text resize option outside of browser/OS features.  Software: No ability to resize text
<b><u>1.4.5 Images of Text</u></b> (Level AA) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>501 (Web)(Software)</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b>  Software: <b>Supports</b>	Electronic Docs: Images of text generally not used unless representing a specific user interface component  Software: No images of text
<b><u>2.4.5 Multiple Ways</u></b> (Level AA) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>501 (Web)(Software) – Does not apply to non-web software</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs) – Does not apply to non-web docs</li> </ul>	Electronic Docs: <b>Partially Supports</b>	Electronic Docs: Multiple navigation components are provided to navigate within sections (header, table of contents), but no overall site map



Criteria	Conformance Level	Remarks and Explanations
<b><u>2.4.6 Headings and Labels</u></b> (Level AA) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b>  Software: <b>Supports</b>	Electronic Docs: Labels and headings are descriptive.  Software: Labels and headings are descriptive.
<b><u>2.4.7 Focus Visible</u></b> (Level AA) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b>  Software: <b>Partially Supports</b>	Electronic Docs: Keyboard focus indicator is visible.  Software: Keyboard focus is generally indicated, though in some cases not easy to determine exact focused element (e.g. not able to determine which table cell in highlighted row is focused)
<b><u>3.1.2 Language of Parts</u></b> (Level AA) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs)</li> </ul>	Electronic Docs: <b>Supports</b>  Software: <b>Does Not Support</b>	Electronic Docs: HTML markup labeled as "en-US"  Software: No API to provide language information.
<b><u>3.2.3 Consistent Navigation</u></b> (Level AA) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software) – Does not apply to non-web software</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs) – Does not apply to non-web docs</li> </ul>	Electronic Docs: <b>Supports</b>	Electronic Docs: Navigation elements are consistent.
<b><u>3.2.4 Consistent Identification</u></b> (Level AA) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software) – Does not apply to non-web software</li> <li>• 504.2 (Authoring Tool)</li> <li>• 602.3 (Support Docs) – Does not apply to non-web docs</li> </ul>	Electronic Docs: <b>Supports</b>	Electronic Docs: Components have consistent identification.
<b><u>3.3.3 Error Suggestion</u></b> (Level AA) Also applies to: Revised Section 508 <ul style="list-style-type: none"> <li>• 501 (Web)(Software)</li> <li>• 504.2 (Authoring Tool)</li> </ul>	Electronic Docs: <b>Supports</b>	Electronic Docs: No user input required  Software: Suggestions to resolve error input not usually provided

Criteria	Conformance Level	Remarks and Explanations
<ul style="list-style-type: none"> <li>602.3 (Support Docs)</li> </ul>	Software: <b>Does Not Support</b>	
<p><b><u>3.3.4 Error Prevention (Legal, Financial, Data)</u></b> (Level AA)</p> <p>Also applies to: Revised Section 508</p> <ul style="list-style-type: none"> <li>501 (Web)(Software)</li> <li>504.2 (Authoring Tool)</li> <li>602.3 (Support Docs)</li> </ul>	<p>Electronic Docs: <b>Supports</b></p> <p>Software: <b>Partially Supports</b></p>	<p>Electronic Docs: Does not modify user-controllable data</p> <p>Software: Data entered by the user is checked for errors on input when possible. Undo/redo functions provided. Modified file warning provided when saving. Once file is saved, software provides no way to reverse submission outside of user-managed backup.</p>

**Table 3: Success Criteria, Level AAA**

Notes: Table 3 has been removed because this product has not been tested for conformance to WCAG Level AAA criteria.

# Revised Section 508 Report

## Chapter 3: [Functional Performance Criteria](#) (FPC)

Criteria	Conformance Level	Remarks and Explanations
302.1 Without Vision	Does Not Support	Software requires visual mode of operation
302.2 With Limited Vision	Partially Supports	Light/dark mode options provided, but no alternative text sizes or other contrast modes provided
302.3 Without Perception of Color	Supports	Alternative labeling (shape, text) are provided for elements with color
302.4 Without Hearing	Supports	No audio content in software
302.5 With Limited Hearing	Supports	No audio content in software
302.6 Without Speech	Supports	No audio input required
302.7 With Limited Manipulation	Partially Supports	No simultaneous actions required, but click targets on some controls (e.g., check boxes)
302.8 With Limited Reach and Strength	Supports	Excessive reach/strength not required outside of normal use of computer
302.9 With Limited Language, Cognitive, and Learning Abilities	Partially Supports	No time-limited functionality. User guide included with software. Minimal/no feedback or guidance in software itself.

## Chapter 5: [Software](#)

Criteria	Conformance Level	Remarks and Explanations
501.1 Scope – Incorporation of WCAG 2.0 AA	See WCAG 2.x section	See information in WCAG 2.x section
<a href="#">502 Interoperability with Assistive Technology</a>	Heading cell – no response required	Heading cell – no response required

Criteria	Conformance Level	Remarks and Explanations
502.2.1 User Control of Accessibility Features	Not Applicable	Not platform software
502.2.2 No Disruption of Accessibility Features	Supports	Does not disrupt platform accessibility features
<b>502.3 Accessibility Services</b>	Heading cell – no response required	Heading cell – no response required
502.3.1 Object Information	Partially Supports	Some user interface elements programmatically determinable but not all
502.3.2 Modification of Object Information	Does Not Support	Not generally supported
502.3.3 Row, Column, and Headers	Partially Supports	Table content can be programmatically determined by screen reader (Windows 10 Narrator)
502.3.4 Values	Does Not Support	Range of allowable values not generally programmatically determinable
502.3.5 Modification of Values	Partially Supports	Some input values (e.g., text or numeric input) may be set by assistive technology
502.3.6 Label Relationships	Partially Supports	Some user interface element labels correctly determined programmatically, some are not or with defects
502.3.7 Hierarchical Relationships	Does Not Support	Hierarchical relationships not explicitly defined or programmatically determinable
502.3.8 Text	Partially Supports	Some text content programmatically available but not all/without defects
502.3.9 Modification of Text	Supports	OS native text inputs used which should be compatible with assistive technology
502.3.10 List of Actions	Does Not Support	Generally not available

Criteria	Conformance Level	Remarks and Explanations
502.3.11 Actions on Objects	Partially Supports	Some native controls compatible with assistive technology
502.3.12 Focus Cursor	Partially Supports	Text insertion point, selection attributes generally exposed for native UI components. Focus tracking available in some places but not without defects
502.3.13 Modification of Focus Cursor	Partially Supports	Compatible in some places
502.3.14 Event Notification	Partially Supports	Some event information available through native components (i.e. window resize)
502.4 Platform Accessibility Features	Not Applicable	Not platform software
<b><u>503 Applications</u></b>	Heading cell – no response required	Heading cell – no response required
503.2 User Preferences	Does Not Support	Application does not honor many underlying operating system display preferences (e.g., text size)
503.3 Alternative User Interfaces	Does Not Support	No alternative user interface provided
<b><i>503.4 User Controls for Captions and Audio Description</i></b>	Heading cell – no response required	Heading cell – no response required
503.4.1 Caption Controls	Not Applicable	No audio/video content
503.4.2 Audio Description Controls	Not Applicable	No audio/video content
<b><u>504 Authoring Tools</u></b>	Heading cell – no response required	Heading cell – no response required
504.2 Content Creation or Editing (if not authoring tool, enter “not applicable”)	Not Applicable	Not authoring tool
504.2.1 Preservation of Information Provided for Accessibility in Format Conversion	Not Applicable	Not authoring tool
504.2.2 PDF Export	Not Applicable	Not authoring tool
504.3 Prompts	Not Applicable	Not authoring tool
504.4 Templates	Not Applicable	Not authoring tool

## Chapter 6: Support Documentation and Services

Criteria	Conformance Level	Remarks and Explanations
<b>601.1 Scope</b>	Heading cell – no response required	Heading cell – no response required
<b><u>602 Support Documentation</u></b>	Heading cell – no response required	Heading cell – no response required
602.2 Accessibility and Compatibility Features	Does Not Support	Software does not contain features marked explicitly as accessibility features and does not document compatibility
602.3 Electronic Support Documentation	See <a href="#">WCAG 2.x</a> section	See information in WCAG 2.x section
602.4 Alternate Formats for Non-Electronic Support Documentation	Supports	Software provides PDF and web-based documentation
<b><u>603 Support Services</u></b>	Heading cell – no response required	Heading cell – no response required
603.2 Information on Accessibility and Compatibility Features	Supports	Contact technical support for information on accessibility and compatibility features.
603.3 Accommodation of Communication Needs	Supports	Email and phone support available: ImmunoassayTS@milliporesigma.com 314-376-3560

## Legal Disclaimer

The information contained herein is based upon our current knowledge. Our company reserves the right to modify/change any information herein, at any time. We disclaim that any use of this Information will be fit for any intended purpose. We expressly disclaim any warranty of outcome, functionality or fitness for purpose and any liability arising from use of this information.

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

Merck, Millipore, Belysa and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.  
© 2024 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

