

My M Safety App

Frequently Asked Questions (FAQs) – January 2021

How do I use the app?

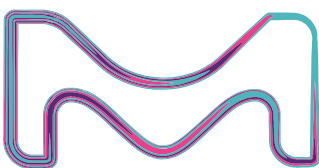
- To get product information, you can use the app's powerful search engine. You can search by name of substance, product or article number, CAS number or use the camera feature in the app to scan the 2D Bar Code on the product label. With just a few additional clicks, your most relevant Safety Data can be looked up or Safety Tags can be printed.

What are the unique features of the app?

- The My M Safety app allows you to easily access Product Safety Information by taking a photo of the 2D Barcode on the product label. Of course, a search by product identifiers is also possible.
- Login options with email or LinkedIn account. This allows you to sync data across multiple devices for easy collaboration.
- Safety Tags containing the most relevant product safety information can be printed directly from the App with a connected printer. This can also be done in bulk mode for multiple products.
- Safety Data is available in multiple languages covering local requirements. The app itself is available internationally with an English and German user interface. We are working on more languages for the user interface as well.
- The app links directly to our Web Shop and customer service. Here, you have access to further product documentation.

What are the newer updates I should look out for? The latest features are...

- Web version available at mysafety.com / mysafety.de
- Apple ID login
- French and Spanish translation of the user interface
- Onboarding / User Guide



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

© 2020 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. Merck and the vibrant M 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. 2020 – 04/05/2020