

Vitroids™ and Lenticule® discs

How to use certified reference materials (CRMs) to perform a spiking protocol for method verification.

Case

Spiking protocol preparation to reach a low level of *Salmonella* inoculation:

3-5 CFU/test portion as requested in ISO 16140-3 (*Protocol for the verification of reference methods and validated alternative methods*) when applying Protocol #3.

Choose an appropriate [CRMs](#), for this case [VT000292 *Salmonella enterica* subsp. *enterica* serovar Abony WDCM 00029](#) is selected, as an example. This product has a mean value range between **15-80 CFU**.

To verify which is the precise value of a specific lot, its [certificate](#) must be checked.

From the certificate of lot BCCG2046, as an example

Sample: <i>Salmonella enterica</i> subsp. <i>enterica</i> serovar Abony WDCM 00029 VT000292		
Certified value (geometric mean value)	Expanded uncertainty (log ₁₀ value)	Expected range
4.9E+01 cfu per disc	0.03	2.4E+01 - 9.8E+01 cfu per disc

In this case the certified value is 49 CFU of *Salmonella* per disc.

Protocol

1. Prepare the initial suspension of the sample to be inoculated (25 g sample + 225 mL diluent).
2. In a separate vial, rehydrate one vitroid in 10 mL Buffered Peptone Water (BPW). Do not spend more than 10-15 minutes for rehydration and spiking (point 3), because the bacteria can start growing in BPW. If the reference material contains 49 cfu per disc (with uncertainty: from 24 cfu to 98 cfu), this *Salmonella* suspension will contain **4,9 cfu/mL** (with uncertainty from 2,4 to 9,8 cfu/mL).
3. Spike the initial suspension (25 g sample + 225 diluent) with 500 microliters of the homogenized *Salmonella* suspension. Considering the reference value, the sample finally would be spiked with a reference value of **2,45 cfu/25 g sample** (with uncertainty between 1,2 and 4,9 cfu/25 g).
It is possible to use a higher amount of *Salmonella* suspension for spiking: for example, if using 800 microliters, the reference value will be still valid (**3,92 cfu/25 g sample**) with better chance to obtain positive results, but the uncertainty of the spiking level will be slightly above the 5 cfu/sample (between 1,92 and 7,84 cfu/25 g).

To place an order or receive technical assistance

Order/Customer Service: **SigmaAldrich.com/order**

Technical Service: **SigmaAldrich.com/techservice**

The Life Science business of Merck operates as MilliporeSigma in the U.S. and Canada.

© 2023 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. Merck, the vibrant M, Millipore, Vitroids and Lenticule 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.

MK_AN12086EN Ver. 1.0 03/2023

Merck KGaA
Frankfurter Str. 250
64293 Darmstadt

