



## Certificate of Analysis

Revvity Health Sciences B.V.  
Rigaweg 22  
9723 TH Groningen  
The Netherlands

Phone: +31-50-544 59 00  
SDS.Groningen@revvity.com  
www.revvity.com

**Product:** Emulsifier-Safe

**Batch #:** 76-23331

**Exp. Date:** March 1, 2026

-----  
The product has been analyzed and is certified to have met the specifications.

Test	Temperature	Results
Appearance		Clear blue fluorescent liquid
Quench parameter		76.0%
Water Holding Capacity	5 °C	≥ 1.1 ml clear
0.01 M PBS Holding Capacity	27 °C	≥ 1.9 ml clear
	20 °C	≥ 2.4 ml clear
Appearance 0.9 ml 1.0 M NaCl	27 °C	Clear within 1.5 minute

Date: August 14, 2023

This Certificate of Analysis has been produced electronically and is valid without a signature. For questions regarding this Certificate of Analysis, please contact your customer service representative.