

Product datasheet

10X Dilution Buffer ab222287

Overview

Product name 10X Dilution Buffer

Product overview 10X Dilution Buffer is available as a stand-alone component for our DCFDA / H₂DCFDA - Cellular ROS Assay Kit ([ab113851](#)).

Prepare 1X Buffer by diluting 10X buffer in ddH₂O.

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	50 ml
10X Dilution Buffer (Sterile)	5 x 10ml

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors