

PRODUCT INFORMATION

SIRT1 FRET Developer

Item No. 10011085



Overview

Contents: The vial contains 50 µl of a lysyl endopeptidase solution.
Storage: -80°C
Stability: ≥2 years

Laboratory Procedures

Thaw the developer on ice. Dilute 15 µl of developer with 735 µl of diluted SIRT1 FRET Assay Buffer (10X) (Item No. 10011080) prior to performing the SIRT1 FRET-Based Screening Assay Kit (Item No. 10010991). The diluted developer is stable for four hours at 4°C.

Precaution

This product may not perform as described if used with other assay reagents or protocols.

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/21/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM