

# PRODUCT INFORMATION



## Resorufin Standard

Item No. 700023

### Overview

---

**Contents:** This vial contains 100  $\mu$ l or 500  $\mu$ l of 2 mM resorufin in DMSO.  
**Storage:** -20°C  
**Stability:**  $\geq$ 2 years

### Laboratory Procedures

---

Follow the applicable kit booklet instructions for preparing the resorufin standard curve.

### Precaution

---

For laboratory research use only. This reagent is not for human or veterinary use.

#### 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, 03/08/2022

#### 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