

PRODUCT INFORMATION



RIPA Buffer Concentrate

Item No. 10010263

Overview

Contents: This vial contains 10 ml of 250 mM Tris-HCl, pH 7.6, containing 750 mM sodium chloride, 5% Tergitol (NP-40), 2.5% sodium deoxycholate, and 0.5% SDS.
Storage: -20°C
Stability: ≥2 years

Laboratory Procedures

Dilute the contents of the vial with 40 ml of HPLC-grade water to yield 50 mM Tris-HCl, pH 7.6, containing 150 mM sodium chloride, 1% Tergitol (NP-40), 0.5% sodium deoxycholate, and 0.1% SDS. This buffer can be used for homogenization of tissue samples or for preparation of cell lysates. The diluted buffer is stable for at least three months 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, 10/11/2018

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