

PRODUCT INFORMATION



Trypan Blue (10X)

Item No. 400292

Overview

Contents: This vial contains 500 µl of Trypan Blue at 0.4% in normal saline.
Storage: Room temperature
Stability: ≥2 years

Laboratory Procedures

To prepare a Quenching Solution, dilute the Trypan Blue 1:10 in Assay Buffer. Add the Quenching Solution to the cells for a couple of minutes and aspirate the excess trypan blue before analysis.

Precaution

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work with the Phagocytosis Assay Kit (IgG FITC) from Cayman (Item No. 500290). 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, 01/11/2022

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