

# PRODUCT INFORMATION

## EDTA Stock Solution (500 mM, pH 8.0)

Item No. 600215



### Overview

---

**CAS Registry No.:** 60-00-4  
**Contents:** This vial contains EDTA (500 mM, pH 8.0).  
**Storage:** 4°C  
**Stability:** ≥1 year

### Laboratory Procedures

---

This bottle contains filtered EDTA (500 mM, pH 8.0). It is ready to use as supplied. For long term storage, we suggest that this product be stored at 4°C; it should be stable for at least one year.

#### 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, 11/12/2025

#### 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](http://WWW.CAYMANCHEM.COM)