

PRODUCT INFORMATION



UltraPure Water

Item No. 400000

Overview

| | |
|-------------------|-------------------------------------|
| Contents: | This vial contains ultrapure water. |
| FW: | 18.0 |
| Storage: | Room temperature |
| Stability: | ≥1 year |

Laboratory Procedures

Cayman's UltraPure water is deionized and free of organic contaminants. It has been purified by reverse osmosis followed by ion exchange and organic scavenging. When bottled, the resistivity of the water was >18.2 MΩ•cm and total organic contaminants were <10 ppb.

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, 01/03/2024

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