

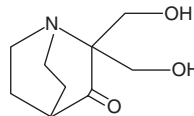
# Product Information



## PRIMA-1

Item No. 63520

**CAS Registry No.:** 5608-24-2  
**Formal Name:** 2,2-bis(hydroxymethyl)-3-quinuclidinone  
**MF:** C<sub>9</sub>H<sub>15</sub>NO<sub>3</sub>  
**FW:** 185.2  
**Purity:** ≥95%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A crystalline solid



For long term storage, we suggest that PRIMA-1 be stored as supplied at -20°C. It will be stable for at least one year.

PRIMA-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the PRIMA-1 in an organic solvent purged with an inert gas. PRIMA-1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PRIMA-1 in these solvents is approximately 25 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of PRIMA-1 can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of PRIMA-1 in PBS (pH 7.2) is approximately 2.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Mutations in the p53 tumor suppressor are found in many different neoplastic cells. PRIMA-1 restores sequence-specific DNA binding and transactivational activity to mutant p53 proteins at relatively high  $\mu$ M to mM concentrations.<sup>1</sup> It is therefore a unique anti-oncogenic substance, working as a re-activator of the apoptotic function of mutant p53 *via* conformational modulation of function-specific epitopes. PRIMA-1 represents a novel lead compound which may be further modified to provide more potent therapeutic agents acting *via* the reactivation mechanism.

### Reference

1. Bykov, V.J.N., Issaeva, N., Shilov, A., *et al.* Restoration of the tumor suppressor function to mutant p53 by a low-molecular-weight compound. *Nature Med.* **8(3)**, 282-288 (2002).

### Related Products

F16 - Item No. 10022 • Pifithrin- $\mu$  - Item No. 10748 • Miltefosine - Item No. 63280 • ICAD (human) Polyclonal Antibody - Item No. 160730 • ICAD (murine) Polyclonal Antibody - Item No. 160735 • CAD (murine) Polyclonal Antibody - Item No. 160740 • Caspase-3 (human) Polyclonal Antibody - Item No. 160745 • Apaf-1 Polyclonal Antibody - Item No. 160780 • Caspase-9 Polyclonal Antibody - Item No. 160790 • Caspase-10 Polyclonal Antibody - Item No. 160795

## Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
custserv@caymanchem.com

**Web**  
www.caymanchem.com

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 07/11/2011