

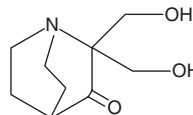
PRODUCT INFORMATION



PRIMA-1

Item No. 63520

CAS Registry No.: 5608-24-2
Formal Name: 2,2-bis(hydroxymethyl)-3-quinuclidinone
MF: C₉H₁₅NO₃
FW: 185.2
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

PRIMA-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the PRIMA-1 in the solvent of choice, which should be purged with an inert gas. PRIMA-1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PRIMA-1 in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

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 solid in aqueous buffers. The solubility of PRIMA-1 in PBS (pH 7.2) is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PRIMA-1 restores sequence-specific DNA binding and transactivational activity to mutant p53 proteins at relatively high μM to mM concentrations.¹ 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. *Nat. Med.* **8(3)**, 282-288 (2002).

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, 11/11/2021

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