

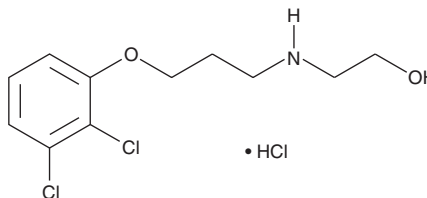
PRODUCT INFORMATION



2,3-DCPE (hydrochloride)

Item No. 10005229

CAS Registry No.: 1009555-55-8
Formal Name: 2-[[3-(2,3-dichlorophenoxy)propyl]amino]ethanol, monohydrochloride
MF: C₁₁H₁₅Cl₂NO₂ • HCl
FW: 300.6
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

2,3-DCPE (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the 2,3-DCPE (hydrochloride) in the solvent of choice, which should be purged with an inert gas. 2,3-DCPE (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2,3-DCPE (hydrochloride) in ethanol is approximately 20 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 2,3-DCPE (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 2,3-DCPE (hydrochloride) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2,3-DCPE is a proapoptotic compound with selectivity for cancer cells versus normal human cells. It induces apoptosis *in vitro* in various cancer cell lines (IC₅₀s = 0.89-2.69 μM) more effectively than in normal human fibroblasts (IC₅₀ = 12.6 μM). 2,3-DCPE down-regulates Bcl-xL, however forced overexpression of Bcl-xL is sufficient to block 2,3-DCPE apoptosis induction.¹ 2,3-DCPE can induce S-phase arrest, upregulate p21, and activate ERK.²

References

1. Wu, S., Zhu, H., Gu, J., *et al.* Induction of apoptosis and down-regulation of Bcl-XL in cancer cells by a novel small molecule, 2[[3-(2,3-dichlorophen-oxy)propyl]amino]ethanol. *Cancer Res.* **64(3)**, 1110-1113 (2004).
2. Zhu, H., Zhang, L., Wu, S., *et al.* Induction of S-phase arrest and p21 overexpression by a small molecule 2[[3-(2,3-dichlorophenoxy)propyl] amino]ethanol in correlation with activation of ERK. *Oncogene* **23(29)**, 4984-4992 (2004).

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, 01/30/2025

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