

DICA

Inhibitor of caspase-3 and -7

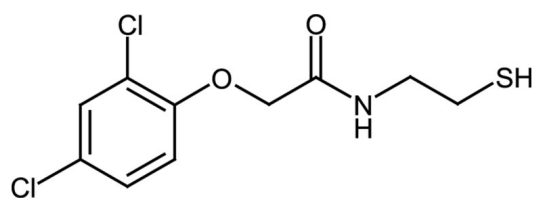
Inhibitor of caspase-3 and caspase-7 binding to an allosteric site and not to the active site.

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Do not store DMSO solutions for more than 1 day.
Handling	Protect from oxygen. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-(2,4-Dichlorophenoxy)-N-(2-mercaptoethyl)acetamide
Appearance	White to off-white solid.
CAS	613665-26-2
Couple Target	Caspase, Caspase-3, Caspase-7
Couple Type	Inhibitor
Formula	$C_{10}H_{11}Cl_2NO_2S$
Identity	Determined by NMR.
MW	280.2
Purity	≥98% (TLC)
Solubility	Soluble in methanol or DMSO.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com