

Cdk/Crk inhibitor

Potent, selective, and ATP-competitive inhibitor of Cdks. Antiproliferative.

Potent, selective, and ATP-competitive inhibitor of Cdks. The Cdk/Crk inhibitor is a broad-spectrum growth inhibitor toward tumor cells.

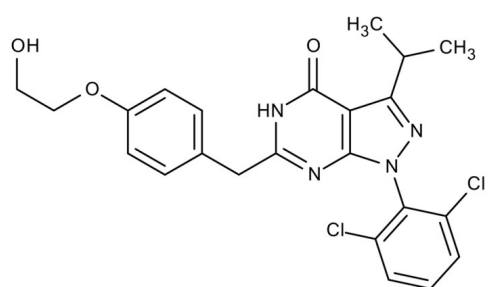
Ordering Information

[Order Online »](#)

ENZ-CHM131-0001	1mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-(2,6-Dichlorophenyl)-1,5-dihydro-6-((4-(2-hydroxyethoxy)phenyl)methyl)-3-(1-methylethyl)-4H-pyrazolo[3,4-d]pyrimidin-4-one, RGB-286147, Cdk7 Inhibitor IV
Appearance	White powder solid.
CAS	784211-09-2
Couple Target	CDK
Couple Type	Inhibitor
Formula	$C_{23}H_{22}Cl_2N_4O_3$
MW	473.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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