

Flavopiridol

CDK inhibitor

Potent inhibitor of cyclin-dependent kinases (CDKs). Induces apoptosis in certain tumor cells. Interacts with multidrug resistance protein 1 (MRP1).

Citations: 26

[View Online »](#)

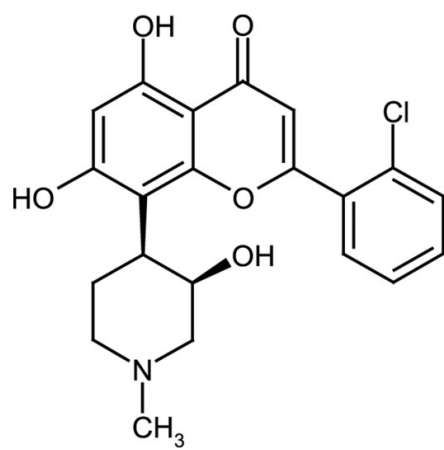
Ordering Information

[Order Online »](#)

ALX-430-161-M005	5mg
ALX-430-161-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Alvocidib, (-)-2-(2-Chlorophenyl)-5,7-dihydroxy-8-[(3R,4S)-3-hydroxy-1-methyl-4-piperidinyl]-4H-1-benzopyran-4-one

Appearance Yellow powder.

CAS 146426-40-6

Couple Target CDK

Couple Type Inhibitor

Formula $C_{21}H_{20}ClNO_5$

Identity Identity determined by 1H-NMR.

MI **14:** 4094

MW 401.8

Purity ≥97% (HPLC)

Solubility Soluble in chloroform; slightly soluble in water.

Source Synthetic.

Last modified: May 29, 2024



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-be@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)