

Diacylglycerol Kinase Inhibitor I (R59 022)

Potentiates PKC activity

Potentiates the activity of protein kinase C (PKC) by inhibiting the diacylglycerol metabolism. Inhibits diacylglycerol kinase without affecting phosphodiesterase or phosphatidyl inositol kinase. It inhibits human platelet DAG kinase in intact cells ($IC_{50}=3.6\mu M$) and in isolated membranes ($IC_{50}=6.7\mu M$).

Citations: 1

[View Online »](#)

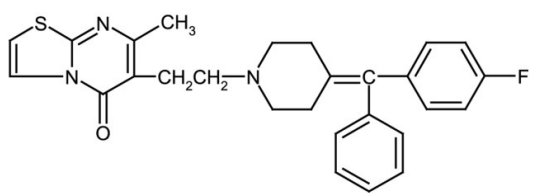
Ordering Information

[Order Online »](#)

ALX-430-027-M005	5mg
------------------	-----

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 6-[2-(4-[(4-Fluorophenyl)phenylmethylene]-1-piperidinyl)ethyl]-7-methyl-5H-thiazolo-[3,2-a]pyrimidine-5-one

Appearance Yellow powder.

CAS 93076-89-2

Couple Target Diacylglycerol kinase

Couple Type Inhibitor

Formula $C_{27}H_{26}FN_3OS$

MW 459.6

Purity ≥99%

Solubility Soluble in DMSO, THF or DMF. Insoluble in water.

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-
be@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)