

BAY 41-2272

Guanylyl cyclase activator

NO-independent activator of soluble guanylyl cyclase (sGC). Activates both isoforms $\alpha_1\beta_1$ and $\alpha_1\beta_2$ of sGC. While the related compound YC-1 (Prod. No. ALX-420-025) also acts as a non-specific phosphodiesterase inhibitor, BAY 41-2272 has no effect on phosphodiesterases. Stimulation of sGC is not blocked by high concentrations of NO scavengers (e.g. PTIO, Prod. No. ALX-430-007) and the combination of BAY 41-2272 with the NO donor DEA NONOate (Prod. No. ALX-430-034) potentiates the activation of sGC. Although BAY 41-2272 alone is not as strong a stimulator of sGC as NO, concentrations as low as 10-100nM stimulate sGC to a level that would be expected to cause biologically important increases in cGMP. ODQ (Prod. No. ALX-270-034), a potent and selective inhibitor of sGC, completely inhibited the effect of BAY 41-2272. The activity of BAY 41-2272 is described in selected literature references.

Citations: 20

[View Online »](#)

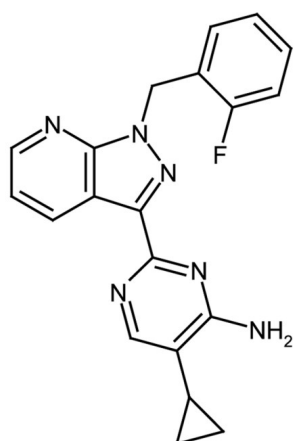
Ordering Information

[Order Online »](#)

ALX-420-030-M005	5mg
ALX-420-030-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 3-(4-Amino-5-cyclopropylpyrimidine-2-yl)-1-(2-fluorobenzyl)-1H-pyrazolo[3,4-b]pyridine

Appearance White to off-white needles.

CAS 256376-24-6

Couple Target Guanylyl cyclase

Couple Type Activator

Formula $C_{20}H_{17}FN_6$

MW 360.4

Purity ≥98% (HPLC)

Solubility Soluble in DMSO, dichloromethane, 100% ethanol or 2-pyrrolidone; practically insoluble in water.



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-
be@enzolifesciences.com](mailto:info-be@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)