

MK-2206 . 2HCl

Highly selective inhibitor of Akt1, Akt2 and Akt3

MK-2206 is a highly selective inhibitor of Akt1, Akt2 and Akt3 with IC₅₀ of 8 nM, 12 nM and 65 nM, respectively.

Citations: 4

[View Online »](#)

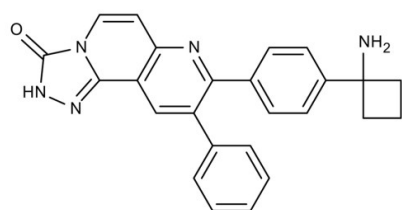
Ordering Information

[Order Online »](#)

ENZ-CHM164-0005	5mg
ENZ-CHM164-0010	10mg

Manuals, SDS & CofA

[View Online »](#)



HCl

HCl

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow powder.
CAS	1032350-13-2
Couple Target	Akt
Couple Type	Inhibitor
Formula	$C_{25}H_{21}N_5O \cdot 2HCl$
Identity	Determined by NMR
MW	480.4
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (14mg/ml) or ethanol (1mg/ml).

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)