

Gardiquimod

TLR7 ligand

Selective ligand for human or mouse Toll-like receptor 7 (TLR7). Induces the activation of NF- κ B in HEK 293 cells expressing TLR7. At high concentrations (3 μ g/ml) slightly activates TLR8. More active than Imiquimod (Prod. No. ALX-420-039).

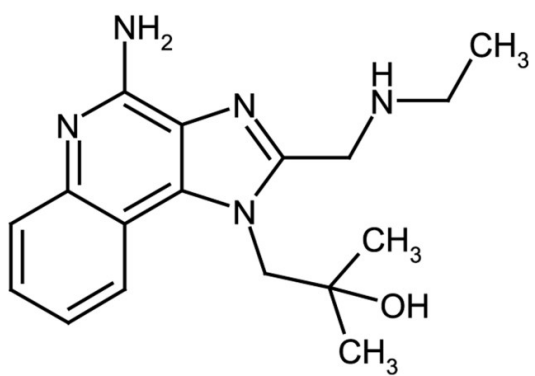
Ordering Information

[Order Online »](#)

ALX-420-040-M025	25mg
ALX-420-040-M100	100mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1-(4-Amino-2-ethylaminomethylimidazo[4,5-c]quinolin-1-yl)-2-methylpropan-2-ol

Appearance White to off-white solid.

Couple Target TLR, TLR7

Couple Type Ligand

Formula $C_{17}H_{23}N_5O$

MW 313.4

Purity ≥98% (HPLC (UV))

Solubility Soluble in DMSO (>40mg/ml), dioxane (7mg/ml), methanol (>40mg/ml) or acetonitrile (2mg/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-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)