

Imiquimod

TLR7 ligand

Topical immune response modifier that inhibits angiogenesis. Up-regulates IL-18 and down-regulates MMP-9 through recognition of Toll-like receptor 7 (TLR7) and subsequent activation of MyD88-dependent pathway.

Citations: 16

[View Online »](#)

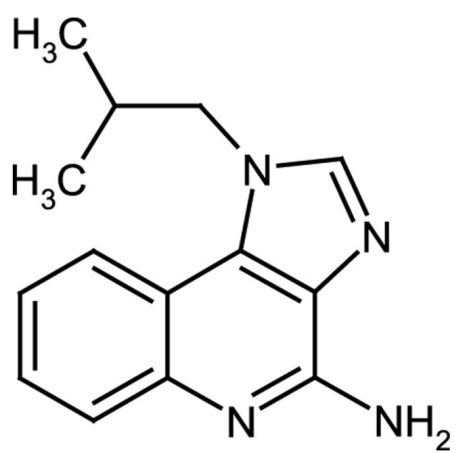
Ordering Information

[Order Online »](#)

ALX-420-039-M100	100mg
ALX-420-039-M250	250mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for up to 1 year after receipt when stored at -20°C. Stock solutions are stable for up to 3 months when stored at -20°C.
Handling	After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-(2-Methylpropyl)-1H-imidazo[4,5-c]quinolin-4-amine, R-837
Appearance	White to off-white powder.
CAS	99011-02-6
Couple Target	TLR, TLR7
Couple Type	Ligand
Formula	$C_{14}H_{16}N_4$
MW	240.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (3mg/ml) or dimethyl formamide; slightly soluble in water (warm to 37°C, vortex); insoluble in 100% ethanol.

Technical Info / Product Notes

Stimulation of TLR7 can be achieved with 1-10µg/ml of the compound for 6 to 24 hours.



ENZO LIFE SCIENCES,
INC.
Phone: +1 949 946 9120
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com