

# 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: 14

[View Online »](#)

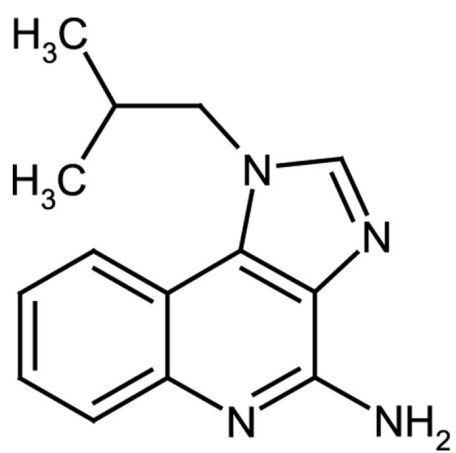
## Ordering Information

[Order Online »](#)

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

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Handling</b>	After reconstitution, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	1-(2-Methylpropyl)-1H-imidazo[4,5-c]quinolin-4-amine, R-837
<b>Appearance</b>	White to off-white powder.
<b>CAS</b>	99011-02-6
<b>Couple Target</b>	TLR, TLR7
<b>Couple Type</b>	Ligand
<b>Formula</b>	$C_{14}H_{16}N_4$
<b>MW</b>	240.3
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in DMSO (3mg/ml) or dimethyl formamide; slightly soluble in water (warm to 37°C, vortex); insoluble in 100% ethanol.
<b>Technical Info / Product Notes</b>	Stimulation of TLR7 can be achieved with 1-10µg/ml of the compound for 6 to 24 hours.



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)