

R-848

Ligand for TLR7 and TLR8

R-848 is a selective ligand for Toll-like receptor 7 (TLR7) in mouse and for TLR7 and TLR8 in human. R-848 displays potent antitumor, proinflammatory, and anti-viral properties. R-848 stimulates antibody secretion, cytokine production (e.g. IL-6, IL-12, and IFN γ), protection from apoptosis, and CD80 upregulation.

Citations: 80

[View Online »](#)

Ordering Information

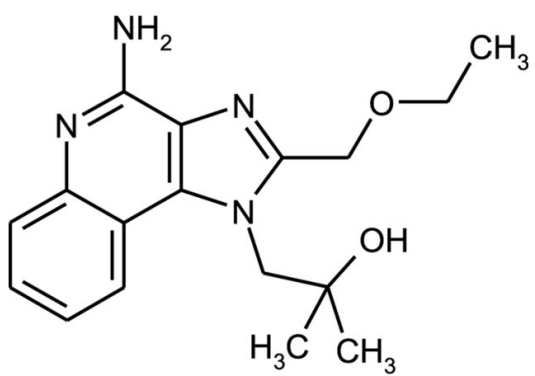
[Order Online »](#)

ALX-420-038-M005	5mg
ALX-420-038-M025	25mg

Manuals, SDS & CoFA

[View Online »](#)

- Highly pure and selective ligand for TLR7 and TLR8
- Highly cited



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 4-Amino-2-(ethoxymethyl)-a,a-dimethyl-1H-imidazo[4,5-c]quinoline-1-ethanol, S 28463, Resiquimod

Appearance White to off-white solid.

CAS 144875-48-9

Couple Target TLR, TLR7, TLR8

Couple Type Ligand

Endotoxin Content Endotoxin-free

Formula $C_{17}H_{22}N_4O_2$

Identity Identity determined by MS, 1H-NMR, and 13C-NMR.

MW 314.4

Purity ≥98% (HPLC)

Solubility Soluble in DMSO, 100% methanol, 5% methanol in dichloromethane; poorly soluble in acetonitrile or ethyl acetate.

Source Synthetic



ENZO LIFE SCIENCES,
Last modified: **May 29, 2024**
Phone: 800.942.0430
info-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-de@enzolifesciences.com

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