

Loxoribine

TLR ligand

Selective ligand for Toll-like receptor 7 (TLR7) via IFN. Activates natural killer cells and primes cytolytic precursor cells for activation by IL-2.

Citations: 8

[View Online »](#)

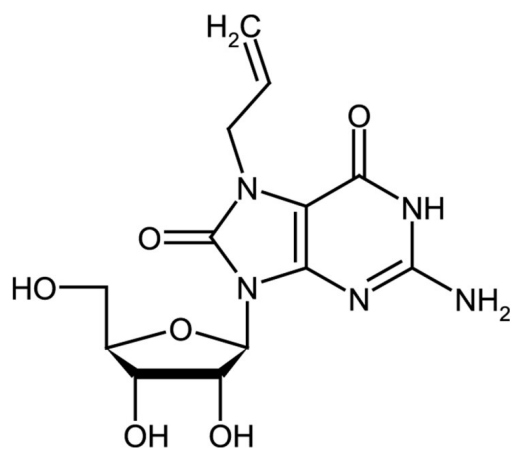
Ordering Information

[Order Online »](#)

ALX-480-097-M025	25mg
------------------	------

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 7-Allyl-8-oxoguanosine, RWJ 21757

Appearance White to off-white solid.

CAS 121288-39-9

Couple Target TLR, TLR7

Couple Type Ligand

Formula $C_{13}H_{17}N_5O_6$

MW 339.3

Purity ≥97% (HPLC (UV))

Solubility Soluble in DMSO (30mg/ml), water (10mg/ml), water:acetonitrile (1:1) (20mg/ml) or water:methanol (1:1) (10mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@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