

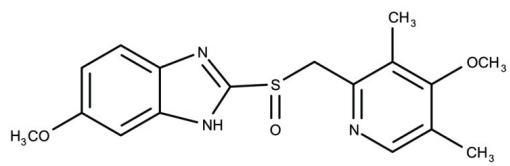
Omeprazole

CYP2C19 inhibitor. H⁺/K⁺ ATPase inhibitor.

Omeprazole inhibits H⁺K⁺-ATPase (IC₅₀ = 5.8 μM). Potent inhibitor of CYP2C19 (K_i=3.1 mM); enhances efficacy of remdesivir (10-fold). Inhibits swelling-dependent chloride channels. Raises stomach pH, which in turn improves oral bioavailability of bioactive peptides such as oxytocin. Displays broad spectrum antiviral activity.

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 2 years after receipt when stored unopened at -20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	73590-58-6
Couple Target	ATPase, Cytochrome P450
Couple Type	Inhibitor
Formula	$C_{17}H_{19}N_3O_3S$
Formulation	Lyophilized.
Identity	Identity determined by NMR.
MW	345.4
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (30 mg/ml) or ethanol (5 mg/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)