

ODN 2216 (TLRGRADE[®]) (synthetic)

TLR9 ligand

CpG oligodeoxynucleotide (Type A) with mixed phosphodiester/phosphorothioate backbone. Specific ligand for human TLR9 (Toll-like receptor 9).

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-746-005-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Aqueous stock solution is stable for 1 day when stored at +4°C.
Handling	Protect from light. For maximum product recovery after thawing, centrifuge the vial before opening the cap. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Couple Target	TLR, TLR9
Couple Type	Ligand
Formulation	Lyophilized. Sterile.
MW	6644 (ammonium salt)
Purity Detail	Activity and endotoxin tested – TLRGRADE®.
Quantity	15.1 nmol
Reconstitution	For a 100µM stock solution, dissolve the total vial content in 151µl endotoxin-free water (included) or PBS (to order separately). To obtain optimal dissolving we recommend the following procedure:- Add 50% of the solvent and let dissolve for 10 min.- Add remaining 50% of the solvent and mix thoroughly.- Moderate warming may aid dissolving.
Sequence	5'-ggGGGACGATCGTCggggggg-3' (lower case letters indicate phosphorothioate linkage).
Source	Synthetic.

Technical Info / Product Notes

Includes 1 vial of ddWater (endotoxin-free) (Prod. No. ALX-505-008).

For the inactive control compound, please see Prod. No. ALX-746-205.

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