

# ODN 2216 (TLRGRADE<sup>®</sup>) (synthetic) (BULK)

## 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-746-055-M001	1mg
------------------	-----

## 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	Activator, Ligand
Endotoxin Content	<0.002 EU/μg (LAL test; BioWhittaker)
Formulation	Lyophilized. Sterile.
MW	6644 (ammonium salt)
Purity Detail	Activity and endotoxin tested – TLRGRADE®.
Quantity	151.1nmol
Reconstitution	For a 200μM stock solution, dissolve the total vial content in 756μl endotoxin-free ddWater (Prod. No. ALX-505-008) or endotoxin-free PBS (Prod. No. ALX-505-007). 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'-ggGGGAC <b>CGATCGT</b> Cgggggg-3'(lower case letters indicate phosphorothioate linkage).
Source	Synthetic.



ENZO LIFE SCIENCES,  
INC.  
Product No. 0430  
info-  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
GmbH AG  
Phone: +41 61 926 8989  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
info-  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

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

Includes 1 vial of ddWater (endotoxin-free) (Prod. No. ALX-505-008).  
For the inactive control compound, please see Prod. No. ALX-746-205.