

# ODN 1585 (TLRGRADE<sup>®</sup>) (synthetic)

## TLR9 ligand

CpG oligodeoxynucleotide (Type A) optimized for NK cell activation with mixed phosphodiester/phosphorothioate backbone. Specific ligand for mouse TLR9 (Toll-like receptor 9).

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Aqueous stock solution is stable for 1 day when stored at +4°C.
<b>Handling</b>	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.
<b>Long Term Storage</b>	+4°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Couple Target</b>	TLR, TLR9
<b>Couple Type</b>	Ligand
<b>Endotoxin Content</b>	<0.002EU/μg (LAL test; BioWhittaker)
<b>Formulation</b>	Lyophilized. Sterile.
<b>MW</b>	6643 (ammonium salt)
<b>Purity Detail</b>	Activity and endotoxin tested – TLRGRADE®.
<b>Quantity</b>	14.9nmol (~99μg). Sufficient for at least 100 cellular activation assays at 0.1-0.5μg/well when performed in a 200μl assay volume (i.e. 96 well microtiter plate).
<b>Reconstitution</b>	For a 100μM stock solution, dissolve the total vial content in 149μl 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.
<b>Sequence</b>	5'-ggGGTCAACGTTGAgggggg-3'(lower case letters indicate phosphorothioate linkage).
<b>Source</b>	Synthetic.

## Technical Info / Product Notes

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

For inactive control compound, please see Prod. No. ALX-746-203.

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)