

# ODN 2243 (TLRGRADE<sup>®</sup>) (synthetic) (Control)

## TLR9 activator

GpC oligodeoxynucleotide with mixed phosphodiester/phosphorothioate backbone.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                  |           |
|------------------|-----------|
| ALX-746-205-T100 | 100 tests |
|------------------|-----------|

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 | <0.002EU/μg (LAL test; BioWhittaker)   |
| Formulation       | Lyophilized. Sterile.  |
| MW                | 6684 (ammonium salt)   |
| Purity Detail     | Activity and endotoxin tested – TLRGRADE®.   |
| Quantity          | 14.8nmol (~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).   |
| Reconstitution    | For a 100μM stock solution, dissolve the total vial content in 148μl 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'-ggGGGAGCATGCTGgggggg-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  
ENZO LIFE SCIENCES  
B.V.  
Phone: +32 20 46 46 46  
info-  
[nl@enzolifesciences.com](mailto:nl@enzolifesciences.com)

France  
ENZO LIFE SCIENCES  
SAS  
Phone: +33 472 440 655  
info-  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
ENZO LIFE SCIENCES  
GmbH  
Phone: +49 7621 5500 526  
info-  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

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

Includes 1 vial of ddWater (endotoxin-free) (Prod. No. ALX-505-008).  
For inactive control compound, please see Prod. No. ALX-746-005.