

# ODN 2395 (TLRGRADE<sup>®</sup>) (synthetic)

## TLR9 ligand

CpG oligodeoxynucleotide (Type C) with phosphorothioate backbone.  
Specific ligand for human and mouse TLR9 (Toll-like receptor 9).

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                  |       |
|------------------|-------|
| ALX-746-020-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   |
| Endotoxin Content              | <0.002 EU/μg (LAL test; BioWhittaker)  |
| Formulation                    | Lyophilized. Sterile.  |
| MW                             | 7068 (ammonium salt)   |
| Purity Detail                  | Activity and endotoxin tested – TLRGRADE®.   |
| Sequence                       | 5'- <b>tcgtcg</b> tttt <b>cggcgcgcgcgcg</b> -3' ( <b>bold</b> letters indicate <b>active</b> , immune-stimulatory portions of CpG sequence, un-bolded letters indicate phosphorothioate linkage) |
| Source                         | Synthetic.   |
| Technical Info / Product Notes | Includes 1 vial of ddWater (endotoxin-free) (Prod. No. ALX-505-008).<br>For the inactive control compound, please see Prod. No. ALX-746-256.   |



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-  
fr@enzolifesciences.com](mailto:info-fr@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)