

ODN 2006 (TLRGRADE[®]) (synthetic)

TLR9 ligand

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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-746-006-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	7720 (ammonium salt)
Purity Detail	Activity and endotoxin tested – TLRGRADE®.
Quantity	12.5nmol (~96µg). 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 125µl 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'- tcgtcgttttgcgttttgcgtt -3'(lower case letters indicate phosphorothioate linkage).
Source	Synthetic.

Technical Info / Product



Includes 1 vial of ddWater (endotoxin-free) (Prod. No. ALX-505-008).
For the inactive control compound, please see Prod. No. ALX-746-206.

ENZO LIFE SCIENCES INC. Phone: 800.942.0430 info-usa@enzolifesciences.com	European Sales Office ENZO LIFE SCIENCES INC. France Phone: +41 61 926 8989 info-eu@enzolifesciences.com	Belgium, France, Germany & Luxembourg Phone: +33 472 440 655 info-be@enzolifesciences.com	Italy Phone: +39 472 440 655 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