

# ODN 1720 (TLRGRADE<sup>®</sup>) (synthetic) (Control)

Negative control for ODN 1668

GpC oligodeoxynucleotide with phosphorothioate backbone.

## Ordering Information

[Order Online »](#)

ALX-746-200-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

Endotoxin Content	<0.002 EU/μg (LAL test; BioWhittaker)
Formulation	Lyophilized. Sterile.
MW	6383 (ammonium salt)
Purity Detail	Activity and endotoxin tested – TLRGRADE®.
Quantity	15nmol (~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 150μl 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'-tccatgagcttcctgatgct-3' (lower case letters indicate phosphorothioate linkage).
Source	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-746-001.



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)