# ODN 2118 (TLRGRADE<sup>®</sup>) (synthetic) (Control)

### **TLR9** activator

GpC oligodeoxynucleotide with mixed phosphodiester/phosphorothioate backbone.

# **Ordering Information**

Order Online »

ALX-746-203-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.

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 6643 (ammonium salt)

Purity Detail Activity and endotoxin tested – TLRGRADE®.

**Reconstitution** 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'-ggGGTCAAGCTTGAggggggg-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-

746-003.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com