

ODN 2118 (TLRGRADE[®]) (synthetic) (Control)

TLR9 activator

GpC oligodeoxynucleotide with mixed phosphodiester/phosphorothioate backbone.

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'-ggGGTCAAGCTTGAgggggg-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.