# ODN 1668 (TLRGRADE<sup>®</sup>) (synthetic)

## **TLR9 ligand**

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

Citations: 5

View Online »

**Ordering Information** 

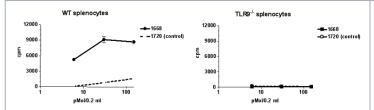
Order Online »

ALX-746-001-C100

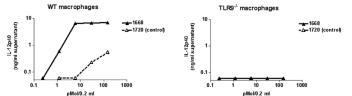
100µg

Manuals, SDS & CofA

**View Online »** 



**Figure 2:** TLR9-dependent mitogenic activity of ODN 1668 (TLRGRADE®) (synthetic) (Prod. No. ALX-746-001) in mouse splenocytes. **Method:** Culture of  $5 \times 10^5$  cells per 200µl well was performed in serum-free conditions after addition of ODNs at the indicated concentrations. Cell proliferation was analyzed after 72 hours with a 16 hours H-thymidine pulse. Only CpG ODN 1668 (Prod. No. ALX-746-001)



**Figure 1:** TLR9-dependent activation of mouse macrophages by ODN 1668 (TLRGRADE®) (synthetic) (Prod. No. ALX-746-001). **Method:** Culture of  $5 \times 10^5$  cells per 200µl well was performed in serum-free conditions after addition of ODNs at the indicated concentrations. Cell activation was measured by IL12-p40 induction. Only CpG ODN 1668 (Prod. No. ALX-746-001)

### **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** <0.002EU/µg (LAL test; BioWhittaker)

**Formulation** Lyophilized. Sterile.

MW 6383 (ammonium salt)

Purity Detail Activity and endotoxin tested – TLRGRADE®.

Quantity 100µg (~15nmol). 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'-tccatgacgttcctgatgct-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-746-200.

Last modified: May 29, 2024

