ODN 2395 (TLRGRADE®) (synthetic)

TLR9 ligand

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

Citations: 1

View Online »

Ordering Information

Order Online »

ALX-746-020-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 7068 (ammonium salt)

Purity Detail Activity and endotoxin tested – TLRGRADE®.

Sequence 5'-tcgtcgttttcggcgcgcgcgcg-3'(bold letters indicate active,

immune-stimulatory portions of CpG sequence, un-bolded

letters indicate phosphorothioate linkage)

Source Synthetic.

Technical Info / Product Notes Includes 1 vial of ddWater (endotoxin-free) (Prod. No.

ALX-505-008).

For the inactive control compound, please see Prod. No.

ALX-746-256.

