Neutral-ODN (Control for iODNs) (endotoxin-free) (synthetic)

Inactive (neutral) control

Inactive (neutral) control ODN for *in vivo* use in rodents (50-150µg per injection) without agonistic nor antagonistic activity.

Several groups have studied the sequence requirements, specificity, signalling pathways and kinetics of the TLR (Toll-like receptor) 9 suppression by 'inhibitory DNA motifs', which led to a revised classification of iODNs, that is independent of the previously thought species preference. Class I: G-stretch ODNs: TLR9-specific competitors, some iODNs may also affect TLR7 and TLR8 signaling

Class II: ODNs with telomeric repeats: TLR-independent inhibitors of STAT signalling (cellular uptake via an "ODN receptor"?)

Class III: Inhibitors of DNA uptake in a sequence independent manner Class IV: Long phosphorothioate ODNs as direct competitors of TLR9 signaling in a sequence independent manner

Slightly modified phosphodiester versions of the most potent inhibitory ODNs were also able to profoundly block the immune activation of macrophages and proved to be valuable tools for in vivo use in experimental animal models of inflammatory and auto-immune diseases.

Citations: 2

View Online »

Ordering Information

Order Online »

ALX-746-256-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

Endotoxin Content <0.002EU/µg

Formulation Lyophilized. Sterile.

MW 4821 (ammonium salt)

Quantity ~20nmol (100μg). Working concentration depends upon

concentration, type and MW of stimulating CpG ODN, the

cellular system and read-out used.

Reconstitution For a 100µM stock solution, dissolve the total vial content

in 270µl 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'-tgctcctggaggggttgt-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).



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