

G-type-iODN (class I) (endotoxin-free) (synthetic)

TLR9 inhibitor

Inhibitory iODN oligodeoxynucleotide with phosphorothioate backbone. Potent sequence of an inhibitory ODN based on poly-G class I sequences targeting TLR9, but **without** additional effect on TLR7 and TLR8 signaling, for *in vivo* use in rodents (50-150µg per injection).

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-746-252-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 | Inhibitor, Ligand |
| Endotoxin Content | <0.002EU/μg |
| Formulation | Lyophilized. Sterile. |
| MW | 6735 (ammonium salt) |
| Quantity | ~15nmol (100μg). Sufficient for ~100 cellular inhibition assays. Working concentration depends upon concentration, type and MW of stimulating CpG ODN, the cellular system and read-out used. Titrate at 3-5 fold molar excess for inhibition. |
| 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'-ctcctattggggttcctat-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 active CpG compounds, please see Prod. No. ALX-746-001.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)