

ODN 1826 (TLRGRADE[®]) (synthetic)

TLR9 ligand

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

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-746-002-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

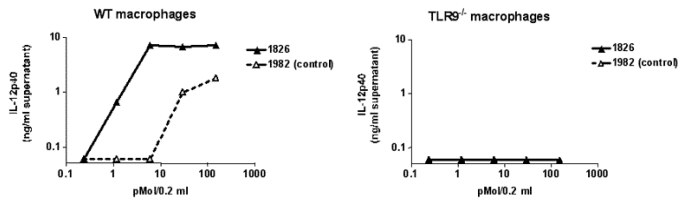


Figure 1: TLR9-dependent activation of mouse macrophages by ODN 1826 (TLRGRADE®) (synthetic) (Prod. No. ALX-746-002). **Method:** Culture of 5×10^5 cells per 200 μ l well was performed in serum-free conditions after addition of ODNs at the indicated concentrations. Cell activation was measured by IL12-p40 induction. Only CpG ODN 1826 (Prod. No. ALX-746-002), but not the corresponding GpC control ODN 1982 (Prod. No. [ALX-746-201](#)) activates wild type mouse macrophages. Activation is absolutely dependent on the presence of TLR9 as addressed with cells derived from TLR9-deficient mice. Note the logarithmic scale for the x and y-axis and the high activity of the active ODNs: max. stimulation is already achieved with 0.010-0.100nmol per 200 μ l well (10-100nM, corresponding to ~0.5-5ug/ml).

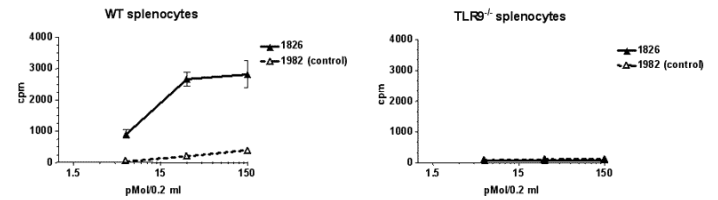


Figure 2: TLR9-dependent mitogenic activity of ODN 1826 (TLRGRADE®) (synthetic) (Prod. No. ALX-746-002) in mouse splenocytes. **Method:** Culture of 5×10^5 cells per 200 μ l well was performed in serum-free conditions after addition of ODNs at the indicated concentrations. Cell proliferation was analyzed after 72 hours with a 16 hours 3 H-thymidine pulse. Only CpG ODN 1826 (Prod. No. ALX-746-002), but not the corresponding GpC control ODN 1982 (Prod. No. [ALX-746-201](#)) induces proliferation of wild type mouse splenocytes. Induction of cell proliferation is absolutely dependent on the presence of TLR9 as addressed with cells derived from TLR9-deficient mice. Note the logarithmic scale for the x and y-axis and the high activity of the active ODNs: max. stimulation is already achieved with 0.010-0.100nmol per 200 μ l well (10-100nM, corresponding to ~0.5-5ug/ml).

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

Alternative Name	CpG-B
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	15nmol (~96μg). 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'-tccatgacggttctctgacggtt-3' (lower case letters indicate phosphorothioate linkage).



Source

Synthetic

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

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

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

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

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

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

Technical Info / Product

Includes 1 vial of ddWater (endotoxin-free) (Prod. No. ALX-505-008).

Notes

For inactive control compound, please see Prod. No. ALX-746-201.

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.