

AT-ODN-1 (endotoxin-free) (synthetic)

TLR9 activator

IFN- γ inducing non-CpG ODN of the AT-type, found in the Malaria genome; TLR9 (Toll-like receptor 9)-dependent immune activation.

CpG ODNs are well known for their immunostimulatory effect mediated by the activation of TLR9. It has been shown that not only CpG ODNs but also some AT-rich DNA fragments (AT-ODNs) are able to exert this effect. AT-ODNs with TLR9-activating and immunostimulatory properties were identified in *Lactobacillus gasseri* as well as in the malarial pathogen *Plasmodium falciparum*. AT-ODNs are useful tools for studying the role of TLR9 in the innate immunological response and in inflammation.

Ordering Information

[Order Online »](#)

ALX-746-024-C100	100 μ g
------------------	-------------

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	Activator, Ligand
Endotoxin Content	<0.0002EU/µg (LAL test; BioWhittaker)
Formulation	Lyophilized. Sterile.
MW	6413
Quantity	15.7nmol
Reconstitution	For a 100µM stock solution, dissolve the total vial content in 157µl 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	TATAATTTTAATTCCAAGA Nucleotides depicted in italics show the corresponding AT-ODN sequence.
Source	Synthetic.
Technical Info / Product Notes	Includes 1 vial of ddWater (endotoxin-free) (Prod. No. ALX-505-008).



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)