

Oligo(deoxyadenosine:deoxythymine) (endotoxin-free) (synthetic)

IFN activator

Strong IFN inducer, **dependent** of IRF3.

Ordering Information

[Order Online »](#)

ALX-746-023-C050	50µg
------------------	------

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

Alternative Name	Oligo(dA:dT)
Couple Target	IFN
Couple Type	Activator
Endotoxin Content	<0.0002EU/μg (LAL test; BioWhittaker)
Formulation	Lyophilized. Sterile.
Quantity	Aliquoted by weight of dry material. Contains polymer, residual salt and water. Sufficient for at least 200 cellular activation assays at 0.01-0.025μg/ml.
Reconstitution	For a 250μg/ml stock solution, dissolve the total vial content in 200μl endotoxin-free buffer or water.
Source	Synthetic.

Technical Info / Product Notes

For *in vitro* stimulation we recommend the following procedure:

1. Bone marrow-derived mouse dendritic cells (BM mDC) are cultured 8×10^4 /100μl in RPMI containing 10% FCS over night at 5% CO₂ and 37°C.
2. The medium is removed and replaced by 100μl OPTI MEM (Invitrogen) and subsequently the transfection mixture is added: 25μl OPTI MEM containing 0.025 to 0.25 Poly(dA:dT) + 25μl OPTI MEM containing 0.5μl Lipofectamine 2000 (Invitrogen) or alternative transfection agents.

3. Over night stimulation at 5% CO₂ and 37°C

4. Measurement of IFN-α/β in culture supernatants



ENZO LIFE SCIENCES
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
(ENZO LIFE SCIENCES)
Phone: +41 61 920 8889
info-eu@enzolifesciences.com

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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