Oligo(deoxyadenosine:de (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.

QuantityAliquoted 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

Last modified: May 29, 2024

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

- 1. Bone marrow-derived mouse dendritic cells (BM mDC) are cultured $8\times10^4/100\mu I$ in RPMI containing 10% FCS over night at 5% CO $_2$ 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.