

Adenovirus type 40 antigen

Native protein.

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-80°C
Shipping	Dry Ice

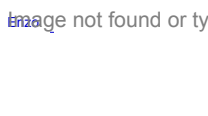
Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Clear liquid.
Formulation	Liquid. In PBS.
Purity	≥99%
Purity Detail	Caesium chloride gradient purified Adenovirus particles
Source	From adenovirus type 40. Inactivated by 2 hours at 54°C.

Last modified: May 29, 2024

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
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