

RecQ DNA helicase (100 µg)

Ordering Information

[Order Online »](#)

CAC-BAM-01-004-EX	100µg
-------------------	-------

Manuals, SDS & CofA

[View Online »](#)

1. Studies on mechanisms of DNA recombination, replication and repair. 2. Application for genetic engineering. For dissociation of special structure DNA.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Solution

Concentration 0.5 mg/ml as measured by BCA method

Formulation 50% glycerol, 20 mM Tris-HCl (pH 7.5), 1 mM EDTA, 50 mM KCl, 1 mM DTT

Purity >90% by SDS-PAGE (CBB staining)

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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