

5X RIPA Buffer with Protease Inhibitor Cocktail

RIPA Buffer extracts proteins from cultured mammalian cells for various downstream applications.

Ordering Information

[Order Online »](#)

CBI-AKR-190	20ml
-------------	------

Manuals, SDS & CofA

[View Online »](#)

- Buffer is able to extract cytoplasmic, membrane, and nuclear proteins
- Suitable for downstream assays such as reporter assays, protein assays, immunoassays and protein purification
- Protease Inhibitor Cocktail included

Handling & Storage

Use/Stability Upon receipt, aliquot and store Protease Inhibitor Cocktail at -20°C and avoid multiple freeze/thaw cycles. Store RIPA Buffer concentrate at room temperature.

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

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