

Brain (rat), (tissue extract)

Brains from male Sprague Dawley rats were suspended in M-PER, Mammalian Protein Extraction Reagent, with protease inhibitor cocktail and homogenized using a Dounce tissue grinder. The tissue lysate was adjusted to 2mg/mL in SDS PAGE sample buffer (40 mM Tris-HCl pH 6.8, 1% SDS, 50 mM DTT, 7.5% glycerol, 0.003% bromophenol blue) and heated for 10 min at 70°C.

Ordering Information

[Order Online »](#)

ADI-LYT-RB100-F	200µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes Western blot control.

Formulation Liquid. In SDS-PAGE sample buffer, pH 6.8, containing 40mM TRIS-HCl, 1% SDS, 50mM DTT, 7.5% glycerol, and 0.003% bromophenol blue.

Source Rat brain extract

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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