

Nuclear extract from HeLa cells

Prepared by high salt extraction of HeLa nuclei (human cervical cancer cell line), this extract is a rich source of HDAC activity, useful for inhibitor screening or immunoprecipitation of HDAC containing complexes. Typically, 0.5µl or less per well is sufficient for assay with the HDAC FLUOR DE LYS[®] fluorometric activity assay (Prod. No. BML-AK500).

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-KI140-0100	100µl
----------------	-------

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

Formulation Liquid. In 0.1M KCl, 20mM HEPES/NaOH, pH 7.9, 20% (v/v) glycerol, 0.2mM EDTA, 0.5mM DTT, 0.5mM PMSF.

Last modified: May 29, 2024

image not found or type not known
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