

HL60 Cell Extract (non-induced)

Immunoblotting standard. Whole-cell extract of human HL-60 leukemia cells. Supplied in SDS-PAGE sample buffer, ready-to-use. Applications: WB control for PARP antibodies, to detect intact 116 kDa PARP. Recommended gel load is 20 µl per lane (~75,000 cells).

Ordering Information

[Order Online »](#)

BML-SW101-0200	200µl
-----------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Formulation Cell extract in SDS-PAGE sample buffer (62.5 mM Tris/HCl, pH 6.8, containing 6 M urea, 10% glycerol, 2% SDS, 0.00125% bromophenol blue and 5% β -mercaptoethanol). The sample is ready to use for SDS-PAGE.



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)