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.

Last modified: May 29, 2024

info-

eu@enzolifesciences.com