

HeLa (heat shocked) (no recovery time), (cell lysate)

Cellular stressors such as heat shock and oxidative stress result in the activation of heat shock protein gene expression by transcriptional regulators such as HSF1. Heat shock proteins play an important role in both normal cellular homeostasis as well as cell survival during and after cellular stress. Heat shock-treated HeLa cells with no recovery time can be a useful tool in studying the regulators of immediate-early gene expression during heat shock stress response. Refer to Ahn and Thiele (2003).

Ordering Information

[Order Online »](#)

ADI-LYC-HL102-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.

Source Human

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)