

# PC-12 (heat shocked) , (cell lysate)

The PC-12 cell line was derived from a transplantable rat adrenal pheochromocytoma. 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.

## Ordering Information

[Order Online »](#)

ADI-LYC-PC101-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.

**Concentration** 2mg/ml

**Formulation** Liquid. In SDS-PAGE sample buffer, pH 6.8, heated for 10 min. at 70°C and containing 40mM TRIS-HCl, 1% SDS, 50mM DTT, 7.5% glycerol, and 0.003% bromophenol blue.

**Source** Rat



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)