

# Hyper Broth™ (single packs)

Hyper Broth™ is an APF Certified™ proprietary media formulation developed by AthenaES™ for improved expression of recombinant proteins in E. coli. The medium is a complex, rich formulation which supplies amino acids, vitamins, and glucose at high levels. This medium yields the highest level of biomass of all AthenaES™ media. The medium is buffered at pH  $7.3 \pm 0.2$  and has the highest buffering capacity of Athena's Expression Media, which permits high glucose levels. For expression of recombinant proteins, induction should be done at cell densities three times that of LB Broth. Cell yields are typically 5 times that of LB Broth depending on the strain. Plasmid copy numbers are significantly higher than they are for LB Broth. This medium is suitable as the base for fermentation cultures and is excellent for feed-batch applications. Biomass yields will depend on the process and the strain employed. Hyper Broth™ powder is shipped with the necessary Glucose Nutrient Mix.

## Ordering Information

[Order Online »](#)

AES-0107-S	10x1L
------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      Ambient

Shipping      Ambient Temperature

**Regulatory Status**      RUO - Research Use Only

## Product Details

Formulation      Powder.



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)