

Power Broth™

Power Broth™ is a proprietary medium 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 a carbon source at higher levels than Terrific Broth⁵. The medium is buffered at pH 7.2±0.2. For expression of recombinant proteins, induction should be done at cell densities three times that of LB Broth. Cell yields are typically 3 to 4 times that of LB Broth in shake flask cultures depending on the characteristics of the strain. The amount of soluble protein accumulated for otherwise insoluble proteins is increased by growing in Power Broth™. This medium is suitable as the base for fermentation cultures. Biomass yields will depend on the process and the strain employed.

Ordering Information

[Order Online »](#)

AES-0106	500g
----------	------

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com