

QuickFold™ Protein Refolding Kit

The AthenaES™ Protein Refolding Kit lets researchers determine the optimal refolding buffer for their proteins. There are many examples of proteins that are recovered from inclusion bodies for both industrial or basic science applications. Methods for purifying inclusion bodies and solubilizing the aggregated protein by denaturation are well established. There is, however, no reliable method for predicting the conditions needed to refold the protein. Thus, the identification of the conditions needed to properly refold the protein remains an empirical science. The purpose of the Protein Refolding Kit is to help simplify the process of identifying the buffer composition and method which is best suited for the refolding of any given protein. Unlike similar products, our refolding kit allows you to examine a wider range of conditions in one experiment.

Ordering Information

[Order Online »](#)

AES-0600	1Kit
----------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Contents 15 Ready-to-Use Refolding Buffers, 10ml each
Dithiothreitol
Glutathione Redox System
Application Manual

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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