

ATP (Energy) Regeneration Solution

This solution contains MgCl_2 , ATP, and ATP regenerating enzymes to recycle ATP hydrolysis products, (AMP, ADP) to ATP. An ideal energy source for conjugation/degradation assays (ubiquitin, proteasome, SUMO, etc.) particularly those that are preformed with crude lysates/impure reagents that may contain ATP-depleting activities.

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-EW9810-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. It is a 10x solution.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Formulation	Liquid. In 50mM TRIS-Cl, pH 7.5.
-------------	----------------------------------

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)