

PRODUCT DATA SHEET



Product name	ATP (Energy) Regeneration Solution (10× solution)				
---------------------	--	--	--	--	--

Product code	EW9810	Batch number	Z05694	Expiry date	12 months from receipt
---------------------	--------	---------------------	--------	--------------------	------------------------

Product information
<p>This solution contains MgCl₂, 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 performed with crude lysates/impure reagents that may contain ATP-depleting activities.</p> <p>Product supplied as 10× solution.</p>

Storage and use
Product is supplied in 50mM Tris-Cl, pH7.5. Store at -80°C.