



Technical Specification Sheet

ver. 1.0 2025/4/17

Full length skeletal myosin II ATPase assay

SKU: CDS086

- Endpoint ATPase format
- High sensitivity
- Rapid turn around

Contacts

Custom Services

tservice@cytoskeleton.com

USA-303-322-2254

Website links

[Myosin general](#)

[Quote form](#)

Parameter	Specification	Notes
Uniprot ID	G1SJQ4	rabbit skeletal muscle MYH3
Amino-acid range	1-1942	
Substrates	F-actin ATP	
Reaction endpoint	ATP → ADP + Pi malachite green + Pi → malachite yellow 650 nm	only endpoint is suitable due to its low activity
Specific activity	80 nmole/mg/min in F-actin with low salt 850 nmole/mg/min in 0.5M KCl	F-actin stimulated ATPase rate. Note full length myosin auto- inhibits its ATPase. 0.5M KCl has been used as a substitute method.
IC50 examples	na	
CV	6%	eight duplicates
Protein source	Cytoskeleton catalog	Cat.# MY02



Additional Information

ATPase raw data

Using the endpoint ATPase assay format, myosin ATPase is accurately measured.

Dose response IC50

4, 8 or 12 pt dose response curves create an accurate IC50 calculation.

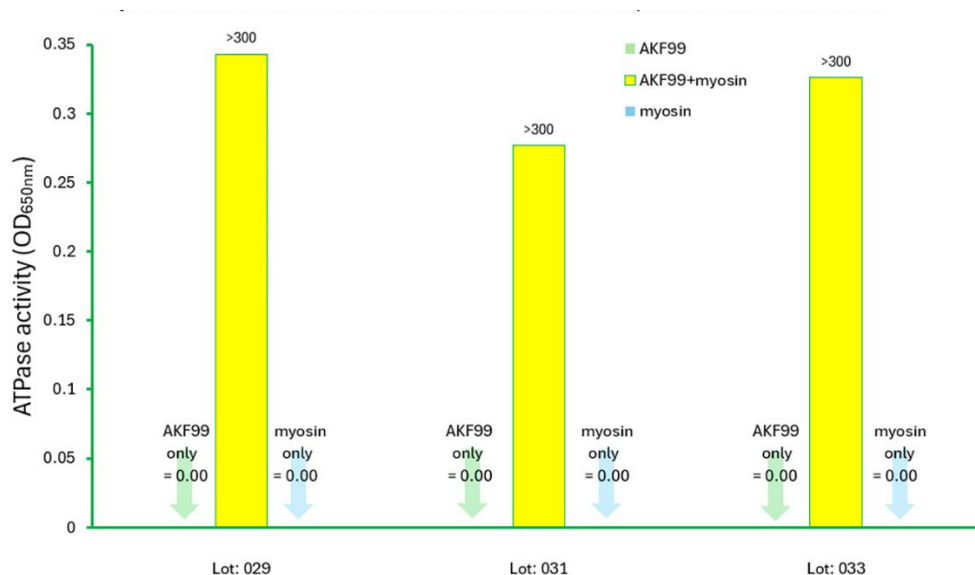
Contact information

Technical Support

Application Support

1830 S Acoma St
Denver, CO 80223.
USA
Phone 303-322-2254
www.cytoskeleton.com

Skeletal myosin ATPase endpoint assay



Legend: Measuring the biological activity of skeletal muscle myosin with pre-formed actin filaments (Cat. #s MY02 and AKF99). The biological activity of myosin ATPase can be determined by an ATPase assay in the presence of actin filaments. Stringent quality control ensures that myosin ATPase is stimulated at least 100-fold over myosin or F-actin alone background control values, as shown in the graph.