

Proteasome 20S

assay kit

A QUANTIZYME® Assay System. Fluorogenic, non-radioactive assay designed to measure chymotrypsin-like protease activity of purified 20S proteasome. This 96-well assay is useful for screening inhibitors and modulators of the 20S proteasome. The kit includes purified, human erythrocyte 20S proteasome (20µg). The Suc-LLVY-AMC fluorogenic peptide substrate supplied with this kit is a proteasome substrate commonly used for measuring chymotrypsin-like peptidase activity. The detection of proteolytic activity is based on the release of free AMC fluorophore and offers the following advantages: non-radioactive, convenient real-time kinetics of substrate cleavage, and excellent sensitivity. A proteasome inhibitor, epoxomicin, is included as a test inhibitor for screening purposes.

Citations: 15

[View Online »](#)

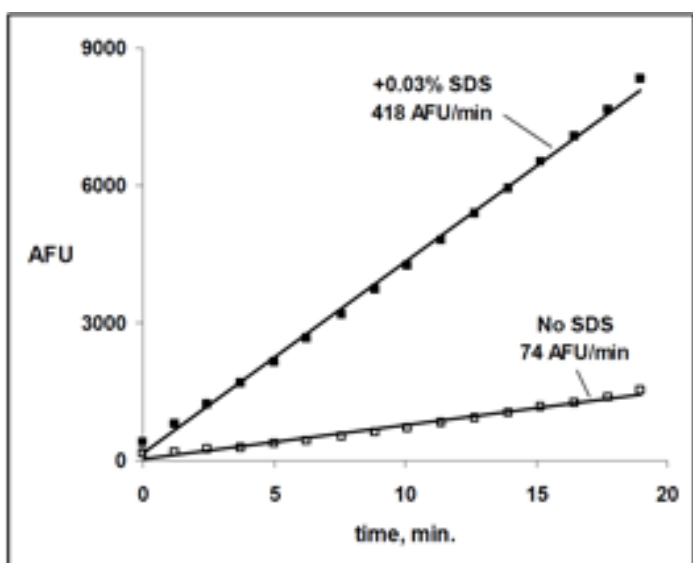
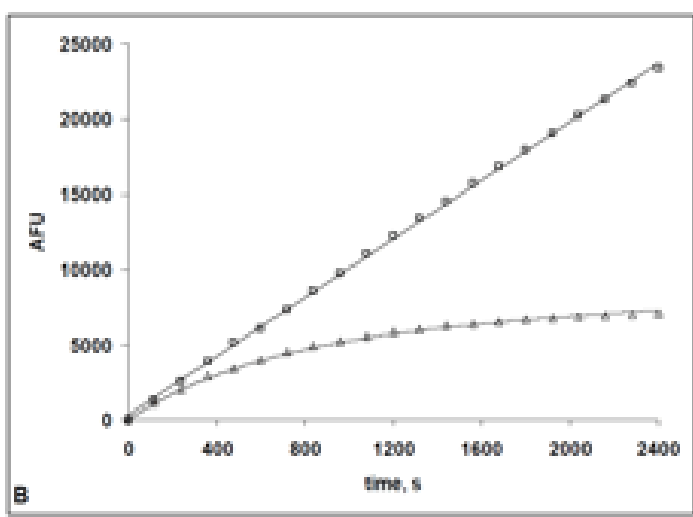
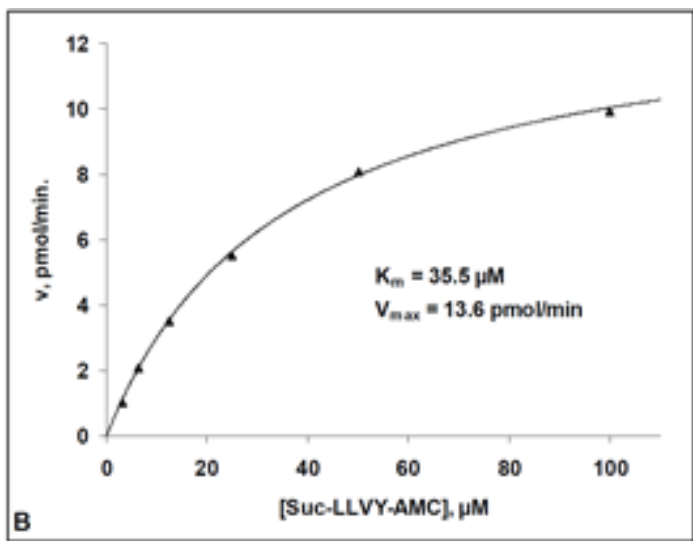
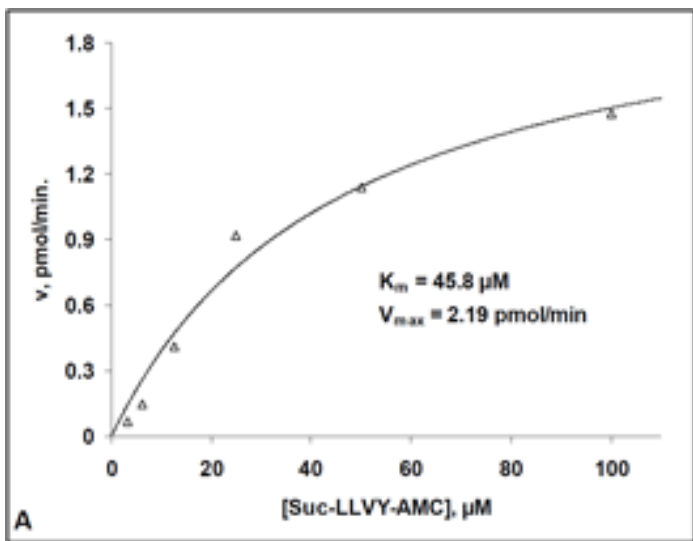
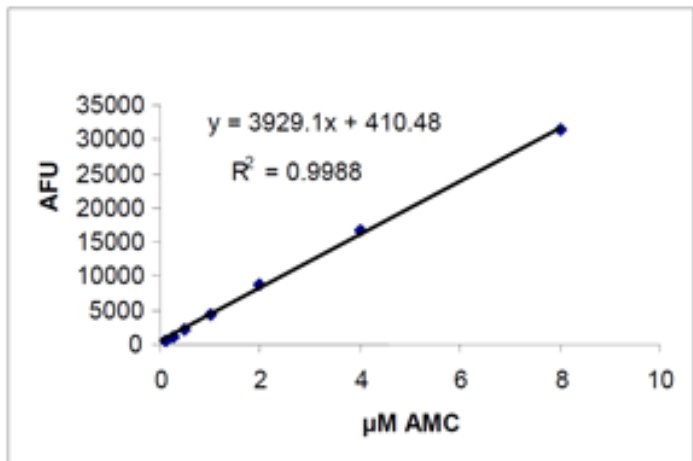
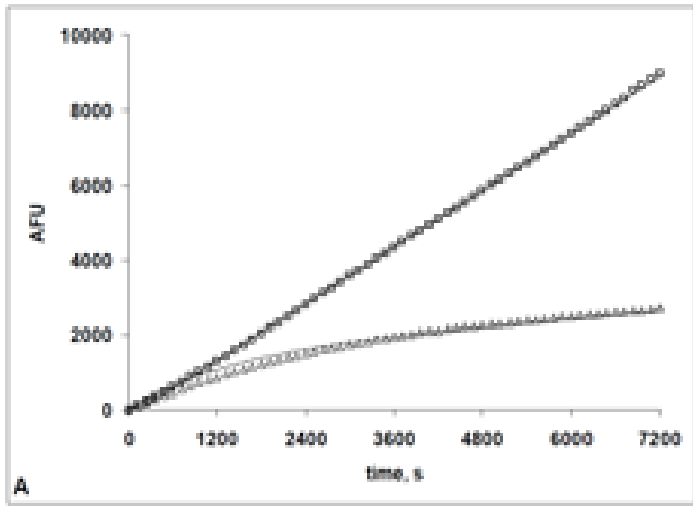
Ordering Information

[Order Online »](#)

BML-AK740-0001	96 wells
----------------	----------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling	Do not store at temperature below recommended storage condition.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	Activity assay, Fluorescent detection, HTS
Contents	Proteasome 20S (human) (purified) (Prod. No. BML-PW8720) Suc-Leu-Leu-Val-Tyr-AMC (Prod. No. BML-P802) Proteasome Assay Buffer (Prod. No. BML-KI340) Sodium Dodecyl Sulfate (Prod. No. BML-KI341) Epoxomicin (Prod. No. BML-PI127) AMC (7-Amino-4-methylcoumarin) Calibration Standard (Prod. No. BML-KI144) 1/2 Volume White Microplates (Prod. No. BML-KI571)

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-
fr@enzolifesciences.com](mailto:info-fr@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)