

BIOMOL® Green

A simple and convenient method for colorimetric phosphate quantitation.

BIOMOL® Green provides a simple and convenient method for colorimetric phosphate quantitation (abs. 600-680 nm). Unlike other molybdate/malachite green-based assays, BIOMOL® Green doesn't require multiple solutions or reagents prepared fresh on the day of the assay. BIOMOL® Green is extremely stable (>6 mos. at 4°C) and is simply mixed, at room temperature, with any enzymatic reaction that has released free phosphate. Reported applications include assays for phospholipid phosphatases, tyrosyl-tRNA synthetase (coupled with pyrophosphatase) and a viral RNA triphosphatase. BIOMOL® Green has perhaps been most widely applied to protein phosphatase assays. Typically, a protein phosphatase (e.g. PTP1B or LAR) is simply incubated with an appropriate phosphopeptide (e.g. IR5,9,10). Addition of BIOMOL® Green stops the reaction and begins color development, which is read 20-30 min. later. BIOMOL® Green may be used in cuvette or microplate-based assays and is ideal for high-throughput applications. Each kit comes with a standardized phosphate solution for assay calibration.

Citations: 159

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-AK111-0250	250ml
BML-AK111-1000	1L

Manuals, SDS & CofA

[View Online »](#)

- Convenient one-step reagent, no mixing needed
- Extremely stable with a long shelf-life, >6 months at +4°C
- Excellent sensitivity, ~50 pmol in 100 µl
- Can be used in cuvette or microplate-based assays, making it ideal for high-throughput applications
- Simple and complete kit that comes with a standardized phosphate solution for assay calibration

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature – Dangerous Good

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, FUNC, HTS

Application Notes Mainly used in protein phosphatase assays. Reported applications include assays for phospholipid phosphatases, tyrosyl-tRNA synthetase (coupled with pyrophosphatase) and a viral RNA triphosphatase.

Contents

Biomol[®] Green Reagent (Prod. No. BML-AK111) (1l or 250ml; liquid in screw cap plastic bottle) Storage: room temperature. Long-term at 4°C; can be stored frozen without deleterious effects.

Phosphate Standard (Prod. No. BML-KI102) (1ml; 800µm phosphate in distilled water) Storage: room temperature. Long-term at 4°C; can be stored frozen without deleterious effects.

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)