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: 156

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 shelflife, >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) (1I 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

