

QuantiFluo™ ALP Reagent (QFALP-6mL)

Liquid Substrate Reagent for Detecting Alkaline Phosphatase Activity

DESCRIPTION

QuantiFluo™ ALP Reagent is a stable 4-methylumbelliferyl phosphate (MUP) liquid substrate reagent formulated for detecting alkaline phosphatase ALP, e.g. in ELISA. This single reagent produces a highly fluorescent product ($\lambda_{\text{ex/em}} = 360/450$ nm) when dephosphorylated by ALP. The fluorescent product can also be visualized under a UV light.

CONTENTS: catalog # QFALP-6mL

QuantiFluo™ ALP Reagent: 6 mL 10x Reagent (pH 10.5)

Instructions for use

The provided reagent is sufficient for 600 assays in 96-well plates. Prior to use, dilute reagent 10-fold in dH₂O. For ELISA, use 100 μ L reagent per well in 96-well plate assays. Incubate 30 min at room temperature. If desired, stop the reaction with 100 μ L phosphate-buffered saline (1x PBS). Read fluorescence at $\lambda_{\text{ex/em}} = 360/450$ nm.

Handling: this product is shipped at ambient temperature. For long-term storage, keep reagent at -20°C. Shelf life of 2 years after receipt.

Bulk orders for volumes >60 mL are available. Please contact info@bioassaysys.com.

LITERATURE

1. Watanabe F, et al (1979). The analysis of alkaline phosphatase isoenzyme using 4-methylumbelliferyl phosphate as substrate on a cellulose acetate membrane. Clin Chim Acta. 91(3):273-6.
2. Gee KR, et al (1999). Fluorogenic substrates based on fluorinated umbelliferones for continuous assays of phosphatases and beta-galactosidases. Anal Biochem. 273(1):41-8.

PRECAUTIONS

Reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents.

Avoid contact and inhalation. Standard laboratory safety procedures should be followed when handling this product. Safety procedures include wearing OSHA approved safety glasses, gloves and protective clothing.

If eye or skin contact occurs, immediately wash affected area with soap and copious amounts of water for at least 15 minutes. Assure adequate flushing by separating the eyelids with fingers. Call a physician to seek medical advice.

If swallowed, wash out mouth with water provided person is conscious and call a physician to seek medical advice.

If inhaled, move individual to fresh air and call a physician to seek medical advice.

TRANSPORT INFORMATION:

DOT (US):	Not regulated
UN Number:	Not regulated
Class:	Not regulated
Packaging Group:	Not regulated

