EvaEZ™ Fluorometric Polymerase Activity Assay Kit

An easy and accurate way to determine the activity of a DNA polymerase without using radioisotopes.



Product attributes

Storage Conditions Store at -10 to -35 °C

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

EvaEZTM Fluorometric Polymerase Activity Assay Kit provides an easy and accurate way to determine the activity of a DNA polymerase without using radioisotopes.

- Ideal for measuring Tag DNA polymerase activity
- Can also be used for other DNA polymerases such as Pfu, Vent, Phusion®, Bst, Phi29, MMLV, AMV, SuperScript®, T4 DNA polymerase, T7 DNA polymerase, Klenow, and E. coli DNA polymerase I
- Kit contains EvaGreen® Dye, a primed template, dNTPs, and MgCl2 in a Tris buffer system
- Available without magnesium by special order

In the presence of DNA polymerase activity, the primer will be extended to form a double stranded product that can bind EvaGreen® Dye, resulting in an increase of fluorescence. The rate of increase of fluorescence is positively correlated to the activity of polymerase. The activity assay can be conducted at temperatures from $4 \,^{\circ}$ C to $75 \,^{\circ}$ C.

EvaEZTM Polymerase Activity Mix is provided at 2X concentration with 5 mM MgCl₂ (final Mg²⁺ concentration in assay is 2.5 mM). Magnesium-free EvaEZTM Polymerase Activity Mix (29051MF) is available on request for customers who wish to control the magnesium concentration in the assay. Contact customer service at order@biotium.com to inquire about availability.

Looking for some advice on using EvaEZ™?

Download our Tech Tip: Protocols and Tips for Success with the EvaEZ™ Fluorometric Polymerase Activity Assay Kit

Phusion and SuperScript are registered trademarks of Thermo Fisher Scientific.

This datasheet was generated on November 6, 2025 at 01:02:32 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/evaeztm-fluorometric-polymerase-activity-assay-kit/