



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## 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

### Product Description

EvaEZ™ 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 Taq 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 MgCl<sub>2</sub> 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 °C to 75 °C.

EvaEZ™ Polymerase Activity Mix is provided at 2X concentration with 5 mM MgCl<sub>2</sub> (final Mg<sup>2+</sup> concentration in assay is 2.5 mM). Magnesium-free EvaEZ™ 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](mailto: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 December 21, 2025 at 03:16:40 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/evaeztm-fluorometric-polymerase-activity-assay-kit/>