

Red hydrogen peroxide assay kit

Red Hydrogen Peroxide Assay Kit provides a simple absorbance-based or fluorometric assay to detect and quantify hydrogen peroxide or peroxidase activity in biochemical assays, cell extracts and in live cells. It can also be used to detect a variety of oxidase activities through enzyme-coupled reactions. The kit is an optimized 'mix and read' assay that is suitable for both 96- and 348-well formats and is compatible with HTS liquid handling.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-51004	1Kit
-----------	------

Manuals, SDS & CofA

[View Online »](#)

- **Flexible:** Quantify hydrogen peroxide in solution, in cell extracts or directly from certain cells
- **Versatile:** Detect a variety of oxidase activities through enzyme-coupled reactions
- **Sensitive:** Detect as little as 10 picomoles of hydrogen peroxide
- **Complete:** Mix and read reagents, with minimal hands-on time
- **Convenient:** Homogenous assay format, fully compatible with HTS automation

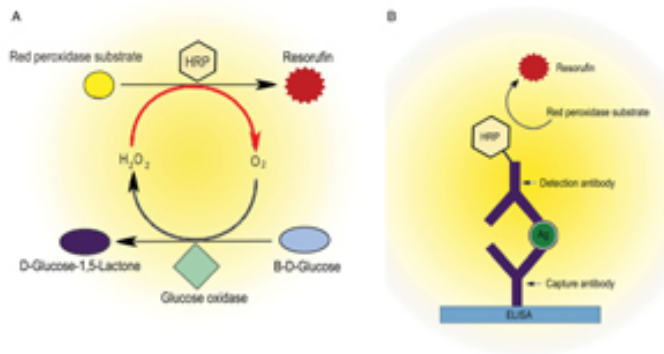


Figure 1: Conversion of Red Peroxidase Substrate into Resorufin. Horseradish peroxidase uses Red Peroxidase substrate as an electron donor during the reduction of hydrogen peroxide to water. The resultant product, Resorufin, is a highly colored and fluorescent compound. (A) Use of assay for detection of glucose oxidase activity. (B) Use of assay in ELISA format to detect HRP-antibody conjugate.

Handling & Storage

Use/Stability Stable for at least one year after receipt when stored as recommended.

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Fluorescence microscopy, HTS

Contents

- Red Peroxidase Substrate** (1 vial)
- Horseradish Peroxidase (HRP)** (20 units)
- Hydrogen Peroxide (H₂O₂) Stock Solution** (200 µl)
- DMSO** (1 ml)
- Assay Buffer** (100 ml)

Quality Control

A sample kit from each lot of Red Hydrogen Peroxide Assay Kit is used to detect H₂O₂ using the protocols described in the user manual. **Minimum Specifications:**

1. A minimum of 10 picomoles of H₂O₂ must be detected in solution under normal conditions.
2. Fluorescence-based detection must yield H₂O₂ specific excitation and emission signals at 570 nm and 585 nm, respectively.
3. Absorbance-based detection must yield an H₂O₂ specific signal at OD₅₇₆.

Quantity For 5 x 96-well plates.

Technical Info / Product Notes

This product is a member of the CELLESTIAL[®] product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL[®] reagents and kits are optimal for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required.