



## Peroxidase Substrate Kit

### Catalog Number 172-1064

**Contents:** Solution A: 180 ml of ABTS color development solution, containing 2,2'-azino-di-[3-ethylbenzthiazoline-6-sulfonic acid].  
Solution B: 20 ml of hydrogen peroxide solution.

**Storage:** The contents of this kit should be stored at 4 °C.

**Shelf life:** 1 year at 4 °C.

- Use:**
1. Allow reagents to come to room temperature before using.
  2. Immediately before use, prepare substrate solution by combining 9 ml of solution A with 1 ml of solution B. Mix well.
  3. After the last rinse following incubation with peroxidase conjugate, add 100  $\mu$ l/well of substrate solution. Gently shake plate during color development to mix contents of wells.
  4. Develop until blue-green color of reaction product appears.
  5. Reaction may be stopped by adding 100  $\mu$ l/well of 2% oxalic acid. Read plate at 415 nm.

This product is intended for research use only. It is not intended for clinical diagnostic purposes.