



Alkaline Phosphatase Substrate Kit

Catalog No. 172-1063

Contents 100 tablets of p-Nitrophenylphosphate, 5 mg each
 100 ml 5x concentrated diethanolamine buffer

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

**Directions
for use**

1. Substrate solution: Immediately before use, mix together 1 ml of 5x diethanolamine buffer and 4 ml of water for each tablet. Make sure each tablet completely dissolves. Substrate solution should be colorless or pale yellow.
2. After last rinse of phosphatase conjugate, add 100 µl well substrate solution. Develop until bright yellow color of reaction product occurs. Read plate at 405-420 nm.
3. Reaction may be stopped by adding 0.4 M NaOH, 100 µl well. Read plate at 405-420 nm.

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