

PRODUCT INFORMATION



HRP Development Pack Item No. 502870

Materials Supplied

Item Number	Item	1 ea Quantity/Size	5 ea Quantity/Size	Storage Temperature
400074	TMB Substrate Solution	1 vial/12 ml	5 vials/12 ml	4°C
10011355	HRP Stop Solution	1 vial/12 ml	5 vials/12 ml	RT

Storage and Stability

Store individual components as directed in the Materials Supplied section and use before the expiration date indicated on the label.

Precautions

TMB Substrate Solution is supplied as a clear solution. If contamination has occurred, the solution may turn blue in color and should not be used. It is sensitive to light and should be stored in the original amber vial.

The HRP Stop Solution in this pack contains an acid. Wear appropriate protection and use caution when handling this solution.

Description

The HRP Development Pack provides a high-sensitivity and low-background chromogenic substrate to measure the enzymatic activity of HRP. When HRP is present, the TMB Substrate Solution reacts with peroxide to form a blue-colored compound that shows peak absorbance at 605 nm. The intensity of this color reflects the level of HRP activity. Adding HRP Stop Solution transforms the blue color to yellow, shifting the absorbance peak to 450 nm. This step halts the reaction and stabilizes the color, allowing precise measurement at 450 nm using a plate reader or spectrophotometer.

ELISA Application

1. Equilibrate the reagents to room temperature before use.
2. Add 100 μ l of TMB Substrate Solution (Item No. 400074) to each well of the plate.
3. Incubate the plate protected from light. Optimal development is obtained using an orbital shaker at room temperature for a minimum of 20 minutes in the dark. *NOTE: The presence of a precipitated product indicates HRP oversaturation and necessitates optimization of the assay conditions, such as enzyme dilution or incubation time.*
4. Add 100 μ l of HRP Stop Solution (Item No. 10011355) to each well of the plate.
5. Wipe the bottom of the plate and read at a wavelength of 450 nm.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/22/2025

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM