

# OPD

34005 34006

0145.2

Number	Description
34005	<b>OPD, 25g</b> <b>Storage:</b> Upon receipt store product at -20°C. Product is shipped at ambient temperature.
34006	<b>OPD Tablets, 50 tablets, 5mg/tablet</b> <b>Storage:</b> Upon receipt store product at 4°C. Product is shipped at ambient temperature.

## Introduction

Thermo Scientific OPD (*o*-phenylenediamine dihydrochloride) is a water-soluble ELISA substrate for horseradish peroxidase. This substrate yields a yellow-orange product with an absorbance maximum at 492nm.

## Procedure for Use

1. Immediately before use, dissolve 5mg of OPD or one tablet in 9mL of water. Add 1mL of Stable Peroxide Buffer (Product No. 34062) to the dissolved OPD. Alternatively, dissolve OPD to 0.5-1.0mg/mL in a buffer containing 0.05M citric acid, 0.05M sodium phosphate; pH 5 and then add 1μL of 30% hydrogen peroxide per 1mL of substrate.
2. Add 100μL of the OPD solution to each microplate well and incubate for 10-30 minutes.
3. For non-stopped assays, measure the absorbance at 450nm.
4. To stop the reaction, add 50-100μL of 2.5M sulfuric acid to each well and measure the absorbance at 490nm.

## Related Thermo Scientific Products

37615	<b>1-Step™ ABTS, 250mL</b>
34024	<b>1-Step Slow TMB-ELISA, 250mL</b>
34022	<b>1-Step Turbo TMB-ELISA, 250mL</b>
34028	<b>1-Step Ultra TMB-ELISA, 250mL</b>
15075	<b>Reagent Reservoirs, 200/pkg</b>
15082	<b>Microtube Racked System, 960 tubes</b>
15036	<b>Sealing Tape for 96-Well Plates, 100/pkg</b>
21124	<b>Streptavidin, Horseradish Peroxidase Conjugated, 2mg</b>
31001	<b>NeutrAvidin® Protein, Horseradish Peroxidase Conjugated, 2mg</b>
37538	<b>StartingBlock (PBS) Blocking Buffer, 1L</b>
37542	<b>StartingBlock (TBS) Blocking Buffer, 1L</b>

---

This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than the original purchaser of the Product ("Buyer").

**No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s).**

There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

Current product instructions are available at [www.thermoscientific.com/pierce](http://www.thermoscientific.com/pierce). For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2011 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.