INSTRUCTIONS

Stable Peroxide Substrate Buffer (10X)

34062

34062

Number Description

Stable Peroxide Substrate Buffer (10X), 100mL

Storage: Upon receipt store product at 4°C. Product is shipped at ambient temperature.

Introduction

The Thermo Scientific Stable Peroxide Substrate Buffer is an ideal formulation for researchers who prefer to make their own horseradish peroxidase (HRP) substrates. The substrate is easily prepared by adding the chosen chromogen to a 1X solution of the Stable Peroxide Substrate Buffer. This buffer is provided at a 10X concentrate, which allows a shelf life of up to one year.

Important Product Information

- A typical concentration of chromogen to use is 1mg/mL; however, specific applications may require a different concentration.
- A precipitate may form when some chromogens are added to the buffer. If a clear substrate solution is desired, filter the solution using a glass-fiber filter, as membrane filters may have affinity for the chromogen.
- Many chromogens, including OPD, ABTS, DAB and TMB dihydrochloride, are readily soluble and can be added directly to the 1X substrate buffer. Some chromogens (e.g., CN, DAB, TMB free base) require dissolving in organic solvents before adding to the 1X substrate buffer.

Substrate Preparation using the Stable Peroxide Substrate Buffer

Add one volume of Stable Peroxide Substrate Buffer (10X) to nine volumes of ultrapure water. Add the chromogen to the 1X-substrate solution immediately before use. Alternatively, dissolve the chromogen in nine volumes of ultrapure water, and then add one volume of the 10X concentrate. Use the substrate immediately after preparation.

Related Thermo Scientific Products

15075 Reagent Reservoirs, 200/pkg 15082 Microtube Racked System, 960 tubes 15036 Sealing Tape for 96-Well Plates, 100/pkg Streptavidin, Horseradish Peroxidase Conjugated, 2mg 21124 NeutrAvidin[®], Horseradish Peroxidase Conjugated, 2mg 31002 StartingBlockTM (PBS) Blocking Buffer, 1L 37538 StartingBlock (TBS) Blocking Buffer, 1L 37542 **EZ-Link[®] Plus Activated Peroxidase Kit** 31489



0407.2



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 <u>www.thermoscientific.com/pierce</u>. 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.