

ABTS

34026

0152.4

Number	Description
34026	ABTS , 50 tablets each containing 10mg of 2,2'-azinobis (3-ethylbenzthiazoline-6-sulfonic acid) diammonium salt with inert filler as the balance CAS #30931-67-0 Storage: Upon receipt store protected from moisture at 4°C. Product is shipped at ambient temperature. Note: ABTS is an irritant causing irritation to eyes, skin, and mucous membranes. Avoid breathing dust. Wear gloves and wash hands thoroughly with water after handling.

Introduction

Thermo Scientific ABTS (2,2'-azinobis [3-ethylbenzthiazoline-6-sulfonic acid] diammonium salt) is a water-soluble ELISA substrate for horseradish peroxidase that yields a green product with absorbance peaks at 410nm and 650nm. Adding sodium dodecyl sulfate (SDS) to the substrate will stop the reaction, but no color change takes place upon stopping.

Procedure for ELISA

A. Reagent Preparation

Note: Stable Peroxide Substrate Buffer (10X) (Product No. 34062), can replace the Substrate Buffer in the procedure; with this buffer, it is not necessary to add peroxide with the ABTS tablet to make the Substrate Solution.

Substrate Buffer	Prepare 100mM phosphate-citrate buffer, pH 5.0 by adding 24.3mL of 0.1M citric acid to 25.7mL of 0.2M dibasic sodium phosphate, then add water to a final volume of 100mL.
Substrate Solution	Add one ABTS tablet to 10mL of Substrate Buffer, then add 10µL of 30% peroxide solution (H ₂ O ₂). Use solution within 1 hour of preparation.
Stop Solution	1% SDS

B. Assay Procedure

1. Add 150µL of Substrate Solution per microplate well.
2. Incubate plate at room temperature for 30-60 minutes.
3. Add 100µL of Stop Solution.
4. Measure absorbance at 405-410nm.

Related Thermo Scientific Products

37615	1-Step™ ABTS, 250mL
34024	1-Step Slow TMB-ELISA, 250mL
34028	1-Step Ultra TMB-ELISA, 250mL

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. 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.