

MATERIALS PROVIDED & STORAGE CONDITIONS

Store the unopened kit at 2-8 °C. Do not use past kit expiration date.

PART	PART #	DESCRIPTION	STORAGE OF OPENED/DILUTED MATERIAL
ELISA Plate-Coating Buffer	896036	60 mL of sterile-filtered 1X PBS for use as a plate-coating buffer in ELISA development kits. Contains no preservatives.	May be stored for up to 3 months at 2-8 °C when handled aseptically.*
Reagent Diluent Concentrate 2 (10X)	841380	2 vials (21 mL/vial) of a 10% BSA solution.	May be stored for up to 1 month at 2-8 °C.* <i>Use and discard diluted Reagent Diluent. Prepare fresh for each assay.</i>
ELISA Stop Solution**	642827	5 vials (12 mL/vial) of Methanesulfonic acid.	May be stored for up to 1 month at 2-8 °C.*
TMB ELISA Substrate	642736	5 vials (12 mL/vial) of substrate.	
Wash Buffer Concentrate	895003	6 vials (21 mL/vial) of a concentrated solution of buffered surfactant with preservative. <i>May turn yellow over time.</i>	
Clear Microplates	640026	5 clear high-binding, flat-bottom polystyrene microplates (360 µL/well). Each microplate consists of twelve removable strips of 8 wells and a frame.	Cover any unused plates with the ELISA Plate Sealers provided.
ELISA Plate Sealers	640197	24 adhesive strips	

* Provided this is within the expiration date of the kit.

** Add 100 µL/well to stop the action of horseradish peroxidase on the substrate, tetramethylbenzidine. If the color in the wells is green or the color change does not appear to be uniform, gently tap the plate to ensure thorough mixing.

REAGENT PREPARATION

Bring all reagents to room temperature before use.

Note: Reagent preparations listed below are sufficient for one 96-well plate. Adjust volumes accordingly.

Reagent Diluent Concentrate 2 (10X):

For assays that require **1% BSA in PBS**, dilute 1:10 in deionized water.

For assays that require **0.1% BSA in TBS**, dilute 1:100 with Tris Buffer + 0.05% Tween® 20, pH 7.2-7.4.

Note: The Reagent Diluent will require additional components for some DuoSets. Refer to the product-specific package insert for more information.

TMB ELISA Substrate - Protect from light.

1X Wash Buffer - If crystals have formed in the concentrate, warm to room temperature and mix gently until the crystals have gone into solution. Add 20 mL of Wash Buffer Concentrate to 480 mL of deionized or distilled water to prepare 500 mL of 1X Wash Buffer.

PRECAUTIONS

The Stop Solution provided with this kit is an acid solution.

Some components in this kit contain a preservative which may cause an allergic skin reaction. Avoid breathing mist.

Wear protective gloves, clothing, eye, and face protection. Wash hands thoroughly after handling. Refer to the SDS on our website prior to use.