

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 3 (5X)	895956	3 vials (21 mL/vial) of a 25% Tween® 20 solution in PBS.	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>	
20X PBS	896034	21 mL of concentrated PBS.	
Clear Microplates	640026	5 clear high-binding, flat-bottom 96-well polystyrene microplates. 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.

1X PBS - Add 3 mL of 20X PBS to 57 mL deionized or distilled water to prepare 60 mL 1X PBS.

Reagent Diluent 3 (diluted 1:5) - Add 12 mL of Reagent Diluent Concentrate 3 to 48 mL of 1X PBS to prepare 60 mL of Reagent Diluent 3 (diluted 1:5) before use.

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 24 mL of Wash Buffer Concentrate to 576 mL of deionized or distilled water to prepare 600 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.