

Wash and Stop Solution for ELISA without Sulfuric Acid

**Code No. MK021 Size: For 96 well plate
x 5 plates**

Description :

This product contains 10X PBS and Tween20 for preparing the wash solution and Stop Solution for stopping the peroxidase reaction, which are used together with TaKaRa Bio's ELISA Kits. Although the Stop Solution is non-sulfuric acid (H₂SO₄) type, it can be used by the same method as 1N H₂SO₄. Use the prepared wash solution and Stop Solution according to the user manual for each ELISA Kit.

Component :

10X PBS	50 ml x 5
Tween20	3 ml
Stop Solution	60 ml

Storage: Room temperature

Preparation :

Wash Buffer

- When PBS is used as wash buffer :

Dilute 50 ml of 10X PBS by adding 450 ml of distilled water to make 1X PBS.

The diluted solution is used as wash buffer after mix well (just before use).

- When 0.1% Tween20/PBS is used as wash buffer :

Dilute 50 ml of 10X PBS by adding 450 ml of distilled water to make 1X PBS.

Add 500 μ l of Tween20 to 500 ml of 1X PBS.

The prepared solution is used as wash buffer after mix well (just before use).

【Note】

Be sure to use the wash buffer specified in the user manual of each ELISA Kit.

Stop Solution

Use directly.

【Note】

As the viscosity of this solution is high , it is necessary to mix well using the plate mixer after adding to the plate.

Precautions for Use :

- * 10X PBS may generate precipitates when stored at low temperature. In case that precipitates are observed, dissolve them completely by warming up in water bath before use.
- * Please be careful not to contact eye and skin with Stop Solution, as it contains weak acid.

Note

This product is for research use only. It is not intended for use in therapeutic or diagnostic procedures for humans or animals. Also, do not use this product as food, cosmetic, or household item, etc. Takara products may not be resold or transferred, modified for resale or transfer, or used to manufacture commercial products without written approval from Takara Bio Inc.

If you require licenses for other use, please contact us by phone at +81 77 565 6973 or from our website at www.takara-bio.com.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product web page. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

All trademarks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions.

