

Version: 01
Update: 03/20/2024

20× Wash Solution

Cat.No. B00063

Description

This Wash Solution is specially formulated and optimized for use in the wash step of ELISA assay.

Product Information

Form	Liquid
Application	ELISA
Recommended Usage	<p>The 20× Wash Solution must be diluted to 1× Wash Solution with deionized or distilled water prior to ELISA wash procedures.</p> <p>Dilute the 20× Wash Solution with deionized or distilled water with a volume ratio of 1:19. For example, dilute 40 mL of 20× Wash Solution with 760 mL of deionized or distilled water to make 800 mL of 1× Wash Solution.</p>
Storage and Stability	<p>The unopened product is stable for at least 12 months from the date of manufacture at 2°C to 8°C, and the opened product is stable for up to 1 month from the date of opening at 2°C to 8°C.</p>
Note	<p>If any precipitate is found in the 20× Wash Solution, incubate the bottle in a water bath (up to 50°C) with occasional mixing until all the precipitate is dissolved.</p>

For research use only. Not intended for human and animal therapeutic or diagnostic use.

生产商：南京金斯瑞生物科技有限公司 江苏省南京市江宁区科学园雍熙路 28 号

Manufacturer: Nanjing GenScript Biotech Co., Ltd. No. 28 Yongxi Road, Jiangning District, Nanjing, Jiangsu, China