

Version: 01
Update: 03/20/2024

Assay Dilution Buffer

Cat.No. B00062

Description

Ready-to-use Assay Dilution Buffer designed to dilute high concentration samples or standards and minimize sample matrix interferences.

Product Information

Form	Liquid
Application	Immunoassay
Recommended Usage	Assay Dilution Buffer can be used directly without dilution. For more information on how to use the Assay Dilution Buffer, please refer to the manuals for the specific kit used.
Storage and Stability	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.
Note	White suspended particles may appear in the reagent, which is normal and can be used.

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