

Version: 01
Update: 03/20/2024

96-well Low Binding White Microplates & Sealers

Cat.No. D00042

Product Information

Form	Microplates and Sealers
Packing unit	5
Plate Color	White
Plate Format	96-well
Plate Feature	12×8-well Strips
Sterile	Non-sterile
Application	Immunoassay
Recommended Usage	For more information on how to use the Microplates and Sealers, please refer to the manuals for the specific kit used.
Storage	Store at room temperature for up to 3 years from date of manufacture. Refer to expiration date on package label.

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