

GenTrack Universal Blockers Quick Manual

(Cat. No.: NGS02406i/ NGS09606i/ NGS38406i)

For Illumina Sequencing Platform

Components	Applicable for NGS Libraries Prepared Using Following Kits*	Storage
GenTrack Universal blockers	GenScript GenTrack Library Preparation Kit	-20°C
	Roche KAPA HyperPrep Kits	
	NEBNext Ultra™ II/IIFS	
	Illumina TruSeq HT/LT DNA Library Prep Kit	

* Please contact us to confirm whether it is applicable for other NGS library preparation kits.

Remain stable for 12 months. Avoid repeated freeze-thaw cycles.

Blocking Guidelines

1. Thaw the Universal Blockers and Human Repeat DNA (HRD, not supplied in this kit) if applicable at room temperature and keep them on ice. Gently vortex and centrifuge briefly.
2. We recommend prepare the DNA Mix in each well or tube as listed below.

Component	DNA Library	HRD*	Universal Blockers [#]
Volume per Reaction	Total 500 ng DNA library	5 µL	2 µL

[#] Recommend to use GenScript GenFisher Hybridization and Wash Kit (Cat. No.: NGS02409/ NGS09609/ NGS38409) together for a better performance.

* HRD is not supplied in the package, and is included in GenFisher Hybridization and Wash Kit.

3. Vortex or pipet the mixture up and down 15~20 times to mix well.
4. Dry down the DNA Mix using a vacuum concentrator (e.g. Eppendorf® Concentrator Plus) set at 45°C, until there is no visible residual liquid.

Welcome to download detailed user guide from GenScript official website.

Version 2019-08-1