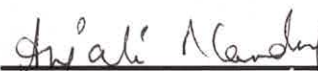


Certificate of Analysis

Product name	Cas9 NLS, S. pyogenes protein with 6x His-tag	Received date	2018/11/06
Cat. No.	GE003 GE004	Detection date	2019/03/10
Lot. No.	GC2CJKFD		
Unit Size	25 µg, 100µg		
Concentration	1 mg/ml		

Quality Control	Specification	Result: Pass/Fail
1. Concentration	1 mg/ml ~ 1.2 mg/ml	Pass
2. Appearance	Clear, colorless	Pass
3. Functional activity	>80% of reference DNA converted to cleavage products	Pass
4. Purity	>90% via SDS-PAGE	Pass
5. Dnase and Rnase check	No detecable Dnase and Rnase activity	Pass
6. Contamination check	No bacterial, fungla, mycoplasma contamination	Pass

Conclusion This lot passes the Quality Control specifications.

Signed by: 
Quality Inspector

Report Date 2019/04/04