

## Certificate of Analysis

### CAT# 381 NanoKAZ-Luciferase protein (NLuc)

Date of Test: 03 April 2018 Lot#E0318

Compound ID: NanoKAZ Luciferase Enzyme  
Quantification: see label, Bradford assay using BSA as standard

Production: recombinant protein produced in E.coli, affinity purified

CAS# Not Registered

Storage: store protein at -80°C,  
avoid repeated freezing

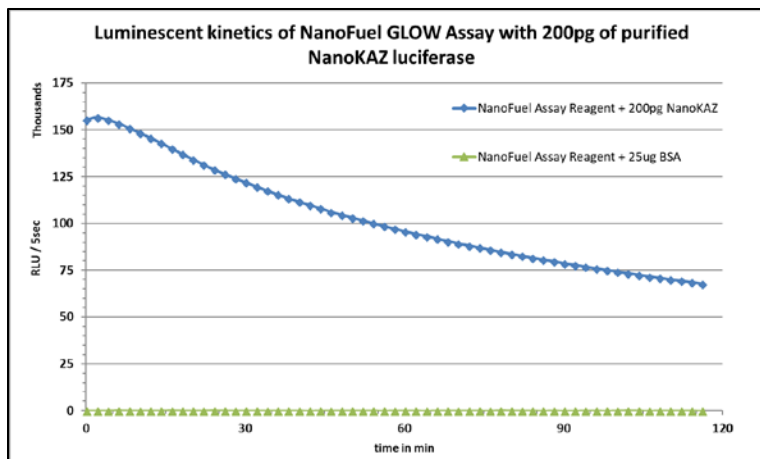
Dilution buffer: NLuc assay buffer from Kit #324 (NLuc FLASH assay) or  
#325 (NLuc GLOW assay), follow instructions in manual

alternative:  
25 mM Tris/Cl, pH 7.8  
150 mM NaCl  
1 mM EDTA  
0.05% (w/v) BSA as stabilizer

Appearance: clear liquid

Activity: recommended CTZ concentration for luminometer: 50uM

NanoKAZ luciferase can be used as a positive control in luminometer assays.



NanoKAZ is patented by US 9,315,783 and US 9,404,145