

CAT# 381 NanoKAZ-Luciferase protein (NLuc)

Date of Test: 9 January 2022 Lot#A1020

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

Production: recombinant protein produced in *E.coli*, affinity purified, stored in TBS, pH 7.8 + 0.1% (w/v) Copolymer

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

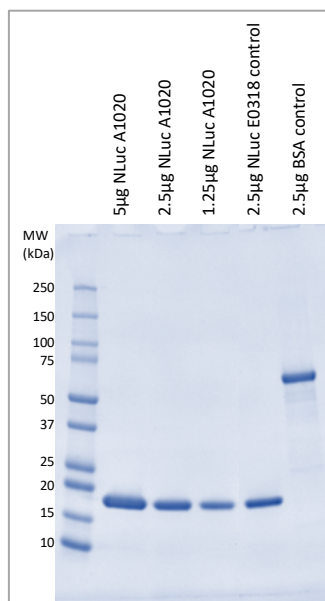
alternative:
TBS or PBS

Appearance: clear liquid

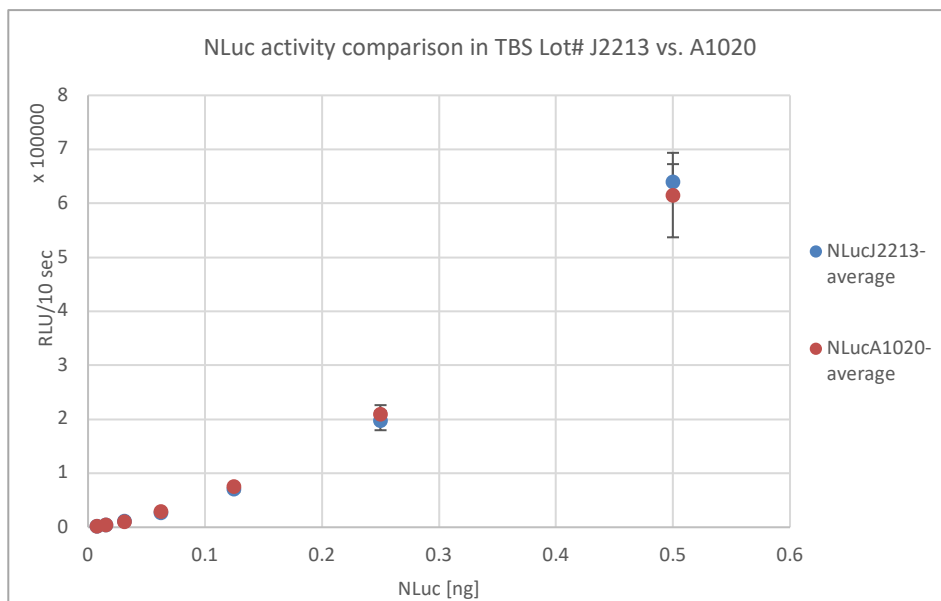
Activity: recommended substrate concentration for luminometer: 50uM

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

SDS-PAGE:



Activity Assay:



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