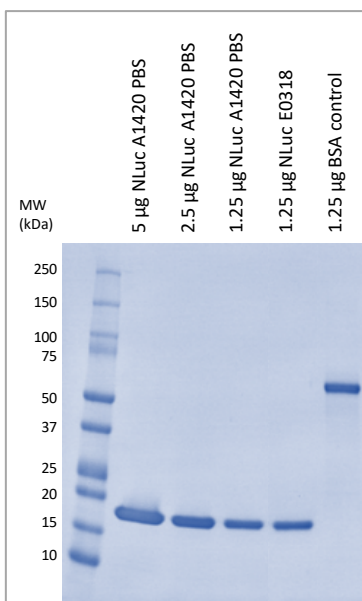


CAT# 3815 NanoKAZ-Luciferase protein (NLuc) in PBS

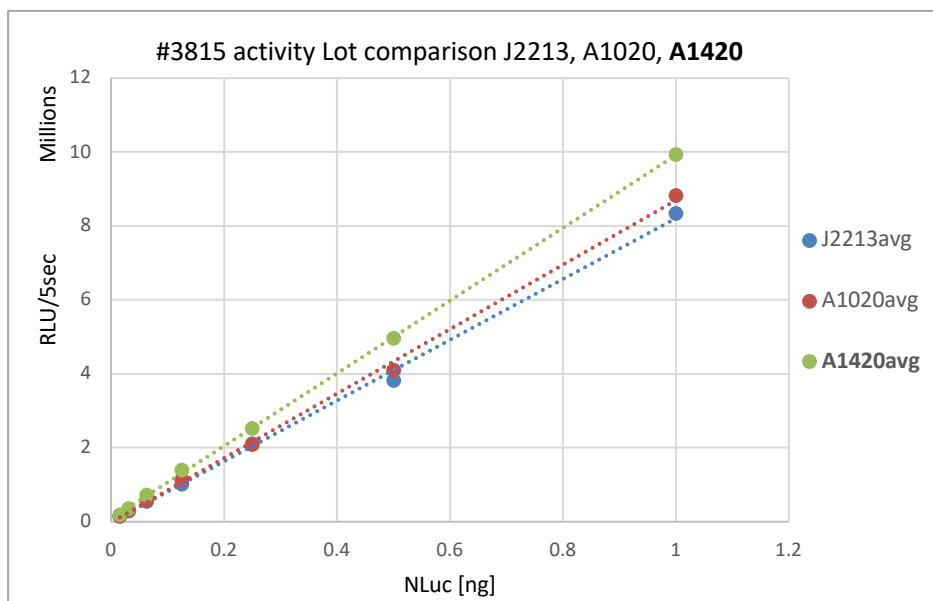
Date of Test:	5 January 2022 Lot#A1020
Compound ID:	NanoKAZ Luciferase Enzyme in PBS
Quantification:	see label, Bradford assay using BSA as standard
Production:	recombinant protein produced in <i>E.coli</i> , affinity purified, stored in PBS, pH 7.4 + 0.1% (w/v) Copolymer
CAS#	Not Registered
Storage:	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 <u>alternative:</u> protein linkage buffer
Appearance:	clear liquid
Activity:	recommended substrate concentration for luminometer: 50uM

NanoKAZ luciferase in PBS can be used for Amine-linked chemistry.

SDS-PAGE:



Activity Assay:



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