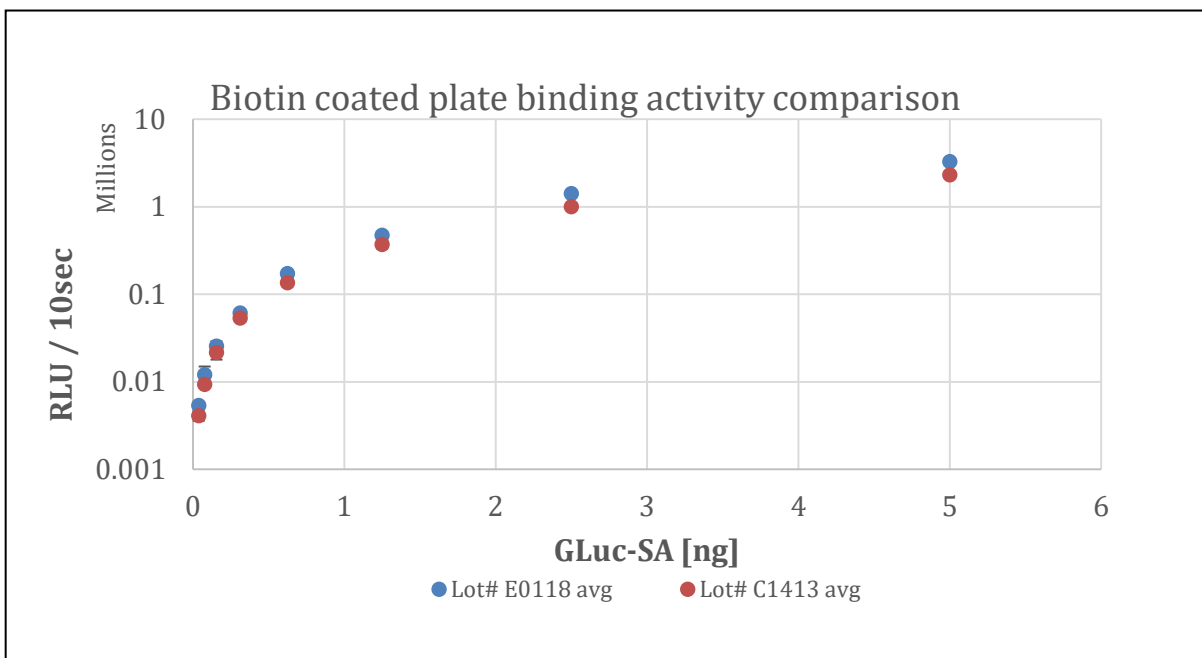


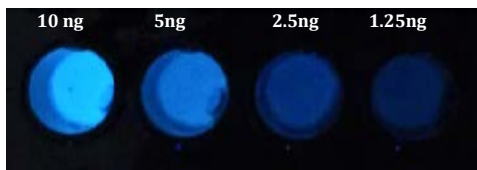
Certificate of Analysis

CAT# 371 Gaussia(M2)-Avitag-biotin-Streptavidin

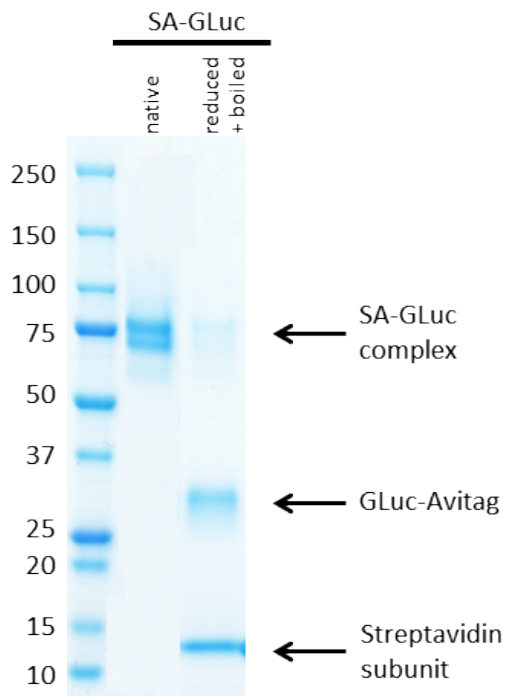
Date of Test:	25 April 2018 Lot#E0118
Compound ID:	Gaussia(M2)-Avitag-biotin-Streptavidin
Quantification:	see label, Bradford assay using BSA as standard
Production:	<i>in vitro</i> fusion of two independently expressed proteins in different organisms
Storage:	store protein at -80°C for maximum shelf-life, avoid repeated freezing
Appearance:	liquid as 1 mg/ml solution in TBS, pH 7.8 + non-proteinogenic stabilizers
Dilution buffer:	TBS or Gaussia dilution buffer (recommended) included in kit Cat.# 319
Activity:	<ul style="list-style-type: none">- recommended CTZ concentration for luminometer: 50µM- around 3-fold brighter than wildtyp GLuc-SA- binding efficiency greater than 90%- very sensitive: use 0.1 to 1 ng of GLuc-SA for biotin coated plates



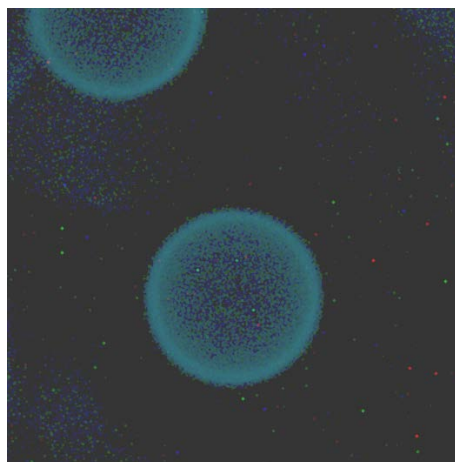
Application examples



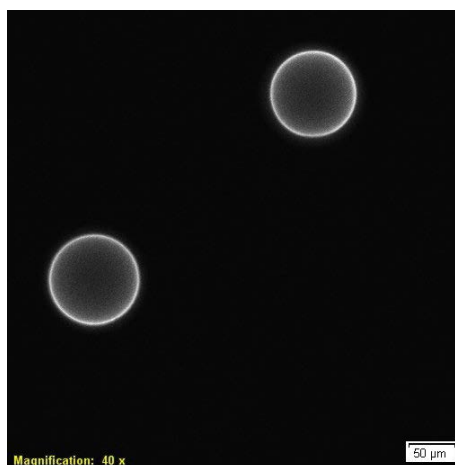
Dilution series (1:2) of GLuc-SA in a black 96-well plate. Picture was taken with a regular camera.



Representative SDS PAGE of GLuc-SA in native and reduced and oxidized form. One GLuc-SA molecule consists of one GLuc-Avitag protein (27 kDa) and 4 Streptavidin subunits (13 kDa).



Biotin-Agarose beads (100 μm) coated with SA-GLuc examined with Nikon microscope with no filter in total darkness.



Biotin-Agarose beads (100 μm) coated with GLuc-SA examined with Olympus LV-200 bioluminescence imaging