

Certificate of Analysis

CAT# 321 Gaussia-Luciferase protein

Date of Test: 06 May 2016 **Lot#G2213**

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

Production: recombinant protein from secretion system lyophilized in proprietary stabilizers

CAS# 481776-15-2
Storage: store dry powder at -20°C or below,
store reconstituted protein at -80°C,
avoid repeated freezing

Reconstitution: as 1 mg/ml solution in TBS, pH 7.8
Dilution buffer: TBS or Gaussia dilution buffer (recommended) included in kit Cat.# 319

Appearance: white lyophilized powder

Activity: recommended CTZ concentration for luminometer: 50uM

SDS Gel Electrophoresis:

1, 5 or 10 ul of 1mg/ml solution of G-Luc was mix with 2x Laemmli buffer (without reducing agent) and directly applied on a 4-20% gel without boiling. Note: reducing the protein will lead to re-oxidation during the SDS-PAGE causing smearing of the protein band if no additional reducing agents are present in the buffer and gel-matrix.

