

Certificate of Analysis

CAT# 321 Gaussia-Luciferase protein

Date of Test: 15 May 2021 Lot#E1221

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: Suggested as 1 mg/ml in TBS, pH 7.8

Dilution buffer: TBS or Gaussia dilution buffer (recommended) included product

Appearance: white lyophilized powder

Activity: tested towards internal standard, recommended CTZ concentration for luminometer: 50-100 μ M

SDS Gel Electrophoresis:

1 μ l of 1mg/ml solution of GLuc was mixed with 2x Laemmli buffer (with and without reducing agent), boiled and directly applied onto a 4-20% SDS gel.

Note: BSA-Fraction V was used as purity control (>98.5%). Molecular weights of a marker standard are listed on the left.

