

# PRODUCT INFORMATION



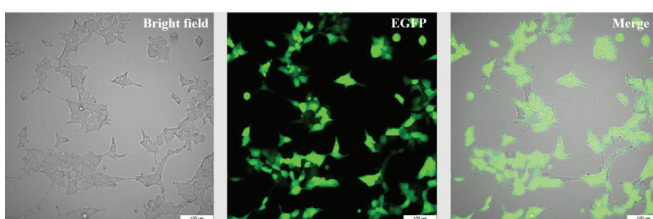
## EGFP mRNA (Cap-1; m1Ψ)

Item No. 45905

### Overview and Properties

<b>Synonym:</b>	Enhanced GFP mRNA
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Supplied as:</b>	A liquid
<b>Storage Buffer:</b>	1 mM sodium citrate, pH 6.4
<b>Concentration:</b>	1 mg/ml
<b>Ex./Em. Max:</b>	489/511 nm

### Image



EGFP mRNA (Cap-1; m1Ψ) 24 hours post transfection in 293T cells.

### Description

EGFP mRNA (Cap-1; m1Ψ) encodes enhanced GFP (EGFP), a variant of *A. Victoria* GFP containing chromophore mutations that increase the sensitivity of the protein and displays excitation/emission maxima of 489/511 nm, respectively.<sup>1</sup> EGFP mRNA (Cap-1; m1Ψ) is polyadenylated and modified with N<sup>1</sup>-methylpseudouridine (m1Ψ) to reduce the host cell immune response and enhance mRNA stability. EGFP mRNA (Cap-1; m1Ψ) can be used for mRNA delivery, cell viability assays, and *in vivo* imaging.

### Reference

1. Zhang, G., Gurtu, V., and Kain, S.R. An enhanced green fluorescent protein allows sensitive detection of gene transfer in mammalian cells. *Biochem. Biophys. Res. Commun.* **227(3)**, 707-711 (1996).

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/22/2026

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM