

PRODUCT INFORMATION

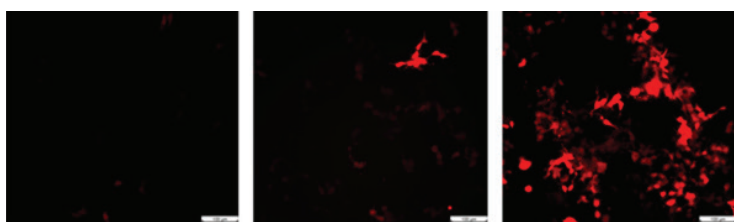
mCherry saRNA (Cap-1; 5mCTP)

Item No. 45906

Overview and Properties

Synonyms:	mCherry Self-amplifying RNA
Storage:	-80°C (as supplied)
Stability:	≥1 year
Supplied as:	A liquid
Storage Buffer:	RMase-free water
Concentration:	1 mg/ml
Ex./Em. Max:	587/610 nm

Image



mCherry saRNA (Cap-1; 5mCTP) fluorescent detection in 293T cells.

Description

mCherry self-amplifying RNA (saRNA) (Cap-1; 5mCTP) encodes mCherry, a red fluorescent protein and derivative of *Discosoma* red fluorescent protein (DsRed) that displays excitation/emission maxima of 587/610 nm, respectively.¹ It is capped with the Cap-1 structure to increase translation efficiency and modified with 5-methylcytidine-5'-triphosphate (5mCTP) to reduce the host cell immune response. mCherry saRNA (Cap-1; 5mCTP) is based on the self-replicating Venezuelan equine encephalitis virus (VEEV), which results in stable expression of mCherry. It can be used for cell labeling *in vitro* and imaging studies in animal models.

Reference

1. Shaner, N.C., Campbell, R.E., Steinbach, P.A., *et al.* Improved monomeric red, orange and yellow fluorescent proteins derived from *Discosoma* sp. red fluorescent protein. *Nat. Biotechnol.* **22**(12), 1567-1572 (2004).

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, 05/20/2026

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