

Transfectamine™ 5000 Transfection Reagent

Catalog number: 60019, 60020, 60021, 60022
Unit size: 50 µL, 0.5 mL, 1 mL, 5 mL

| Component | Storage | Amount (Cat No. 60019) | Amount (Cat No. 60020) | Amount (Cat No. 60021) | Amount (Cat No. 60022) |
|---|--|------------------------|------------------------|------------------------|------------------------|
| Transfectamine™ 5000 Transfection Reagent | Freeze (< -15 °C), Minimize light exposure | 1 vial (50 µL) | 1 vial (0.5 mL) | 1 vial (1 mL) | 1 bottle (5 mL) |

OVERVIEW

Transfectamine™ 5000 Transfection Reagent is a powerful and versatile transfection reagent for the introduction of nucleic acids into eukaryotic cells, or more specifically, into animal cells. It can effectively transfect a variety of payloads into a variety of adherent and suspension cell lines. It can be used for plasmid DNA transfection as well as siRNA- and shRNA-based gene knockdown experiments and gene expression studies. It offers consistently high transfection efficiency in a wide variety of adherent and suspension cell lines, including difficult-to-transfect cells. The low toxicity of Transfectamine™ 5000 also allowed higher viability of transfected cells. Transfectamine™ 5000 is easier to use compare to most other transfection reagents and does not require special medium.

AT A GLANCE

Protocol Summary

1. Prepare cells for transfection
2. Prepare Transfectamine™ 5000-DNA mixture
3. Add Transfectamine™ 5000-DNA mixture to cell culture
4. Culture overnight
5. Analyze transfection efficiency with appropriate method

Important Note

Thaw component at room temperature before starting the experiment.

PREPARATION OF WORKING SOLUTION

1. Mix 2.5 µg of DNA with 200 µL of serum-free medium
2. Add 7.5 µL of Transfectamine™ 5000 to Step 1
3. Mix well and incubate at room temperature for 20 minutes.

Note: Ratio of Transfectamine™ 5000 and DNA need to be optimized for different cell line, in general: Transfectamine™ 5000 Transfection Reagent (µL) to DNA (µg) Ratio = 3 - 5 µL to 1 µg

Sample protocol detail for 6-well and 10 cm plate

| Component | 6 well plate (per well) | 10 cm plate |
|---|-------------------------|-------------|
| Fresh culture medium | 2 mL | 6 mL |
| Plasmid | ~2.5 µg | ~7.5-10 µg |
| Serum-free medium | 200 µL | 600 µL |
| Transfectamine™ 5000 Transfection Reagent | ~7.5 µL | ~22.5 µL |

SAMPLE EXPERIMENTAL PROTOCOL

Preparation of Cell Culture

1. Culture cells to ~ 90% confluency at time of transfection.
2. Replace with fresh growth medium before transfection. For example, replace with 2 mL of medium per well for 6-well plates and 6 mL of medium for 10 cm plates.

Transtection Protocol

1. Add Transfectamine™ 5000 -DNA mixture to culture plate and culture overnight.

Note: Recombinant protein can start to be detected as early as 16 hours post-transfection. Maximal expression level may be observed 72~96 hours post-transfection.

EXAMPLE DATA ANALYSIS AND FIGURES

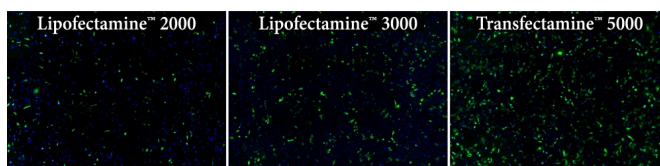


Figure 1. Transfection efficiency comparison in HeLa cells using Transfectamine™ 5000, Lipofectamine 2000 and Lipofectamine 3000 reagents. Each reagent was used to transfect HeLa cells in a 96-well format, and GFP expression was analyzed 24 hours post-transfection. Transfectamine™ 5000 transfection reagent provided higher GFP transfection efficiency compared to Lipofectamine 2000 and Lipofectamine 3000 reagents.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.