

## User Manual

# Alkaline Phosphatase Detection Kit

For 100 tests

**SCR004****FOR RESEARCH USE ONLY****Not for use in diagnostic procedures. Not for Human or Animal Consumption.**

## Product Overview

Embryonic stem (ES) cells are totipotent cells derived from the inner cell mass (ICM) of preimplantation mammalian embryos and are capable of unlimited, undifferentiated proliferation *in vitro*.<sup>1,2,3</sup> Undifferentiated murine ES cells can be maintained *in vitro* for extended periods in media containing the cytokine, leukemia-inhibitory factor (LIF) or our proprietary ES cell culture reagent, ESGRO®.<sup>4,5</sup> The undifferentiated state of ES cells can be characterized by high level of expression of Alkaline Phosphatase (AP),<sup>6</sup> the expression of surface markers including SSEA and TRA antigens and the transcription factor Oct-4.

Our Alkaline Phosphatase Detection Kit (SCR004) is a specific and sensitive tool for the phenotypic assessment of ES cell differentiation by the determination of AP activity.

Available separately from us are the monoclonal antibodies TRA-2-49 (MAB4349) and TRA-2-54 (MAB4354), which permit the detection of Liver/Bone/Kidney isozyme of AP as expressed by ES cells. When used in flow cytometry, these reagents permit a quantitative assessment of AP expression by ES cells.<sup>7</sup>

## Materials Provided

- Fast Red Violet solution (0.8 g/L stock) (90239). Two 15 mL bottles.
- Naphthol AS-BI phosphate solution (2 mg/mL) in a buffer pH 8.5 (CS235583). Two 15 mL bottles.

## Materials Required (Not supplied)

- Fixative (e.g., 4% Paraformaldehyde)
- 1X Rinse Buffer (e.g., TBST: 20 mM Tris-HCl, pH 7.4, 0.15 M NaCl, 0.05% Tween®-20)
- Hematoxylin (optional)
- Microscope

## Storage and Stability

The Alkaline Phosphatase Detection Kit consists of two components used for determining AP activity.

When stored at 2-8 °C, the kit components are stable up to the expiration date. Do not freeze or expose to elevated temperatures. Discard any remaining reagents after the expiration date.

## Protocol

### Preparation of Reagents

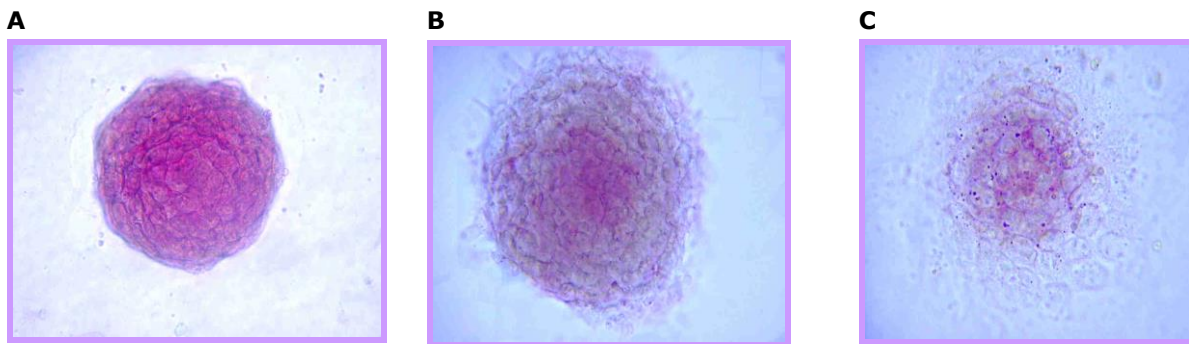
**Naphthol/Fast Red Violet Solution:** Mix Fast Red Violet (FRV) with Naphthol AS-BI phosphate solution in a 1:1 ratio (FRV:Naphthol).

### Staining Protocol

#### Alkaline Phosphatase Staining Procedure

1. Culture ES cells for five days prior to analyzing AP activity, at low to medium density.  
**Note:** This time-period is critical if activity levels of AP needs to be observed. According to our findings, five days of culturing are optimal for good AP stain visualization.
2. On day five, aspirate media and fix ES or EC cells with a fixative (e.g., 4% Paraformaldehyde in PBS) for 1-2 minutes.  
**Note:** Do not over fix. Fixing cells longer than 2 minutes will result in the inactivation of alkaline phosphatase.
3. Aspirate fixative and rinse with 1X Rinse Buffer. DO NOT allow wells to dry.
4. Prepare reagents for Alkaline Phosphatase staining as described in "[Preparation of Reagents](#)" section.
5. Add enough stain solution to cover each well (e.g., 0.5 mL for a well of a 24-well plate). Incubate in dark at room temperature for 15 minutes.
6. Aspirate staining solution and rinse wells with 1 X Rinse Buffer. Cover cells with 1 X PBS to prevent drying and then count the number of colonies expressing AP (red stem cell colonies), versus the number of differentiated colonies (colorless).  
**AP staining criteria:** Greater than 90% of colonies should remain undifferentiated and express alkaline phosphatase in the well containing  $10^3$  Units of LIF. P value shall be  $\geq 0.05$ .

#### Staining with Alkaline Phosphatase Detection Kit



**Figure 1:** Alkaline Phosphatase staining of ES cells: **(A)** High magnification revealed undifferentiated ES cells (mouse MBL.5 cell line)-cultured for five days in media containing our LIF/ESGRO®. A concentration of  $10^3$  Units/mL is used for inhibition of differentiation. **(B)** Differentiated ES cells-cultured at low-medium density for three days in media without any LIF/ESGRO®. **(C)** Differentiated ES cells-cultured at low-medium density for six days in media without any LIF/ESGRO®.

## Related Products

The following related products are available from us as separate items:

- ES Cell Characterization Kit (SCR001)
- ES Marker Sample Kit (SCR002)
- ES Cell 3D Collagen Culture Kit (SCR003)
- SSEA-1 Monoclonal Antibody, purified 100 mg (MAB4301)
- SSEA-3 Monoclonal Antibody, purified 100 mg (MAB4303)
- SSEA-4 Monoclonal Antibody, purified 100 mg (MAB4304)
- TRA-1-60 Monoclonal Antibody, purified 100 mg (MAB4360)
- TRA-1-81 Monoclonal Antibody, purified 100 mg (MAB4381)
- TRA-2-49 Monoclonal Antibody, purified 100 µg (MAB4349)
- TRA-2-54 Monoclonal Antibody, purified 100 µg (MAB4354)
- Murine LIF, 5 mg (LIF2005), 10 mg (LIF2010)
- ESGRO®, 1 x 10<sup>6</sup> units (ESG1106), 1 x 10<sup>7</sup> units (ESG1107)

## References

1. Evans M, Kaufman M. *Nature* 292, 154 (1981).
2. Martin G. *Proc Natl Acad Sci USA* 78, 7634 (1981).
3. Morrison SJ, Shah NM, Anderson DJ. *Cell* 88, 287-298 (1997).
4. Smith AG, Heath JK et al. *Nature* 336, 688-690 (1988).
5. Williams RL, Hilton DJ et al. *Nature* 366, 684-687 (1988).
6. Pease S, Braghetta P et al. *Develop Biol* 141, 344 (1990).
7. Draper J, et al. *J.Anat.* 200, 249 258 (2002).

---

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

### Technical Assistance

Visit the tech service page at [SigmaAldrich.com/techservice](https://SigmaAldrich.com/techservice).

### Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/terms](https://SigmaAldrich.com/terms).

### Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://SigmaAldrich.com/offices).

The life science business of Merck KGaA, Darmstadt, Germany operates as MilliporeSigma in the U.S. and Canada.

MilliporeSigma, ESGRO and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2009-2025 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

Document Template 20769660 Ver 4.0

SCR004 Ver 6.0, Rev 15MAY2025, AV

**Millipore  
Sigma**