

PRODUCT INFORMATION

RHB-Basal™

Catalog Number: Y40000

Size: 500 mL

Applications: Demonstrated applications of RHB-Basal (with addition of supplements) include:

- Maintenance and expansion of adherent mouse and human neural stem (NS) cells
- Neural differentiation of mouse embryonic stem (ES) cells in monolayer culture
- Derivation of NS cells from mouse ES cells and fetal/adult tissues
- Differentiation of mouse and human NS cells into functional neurons
- Maintenance and expansion of glioma neural stem (GNS) cell lines

Description: RHB-Basal is a proprietary, defined, serum-free basal medium specifically formulated for the propagation and differentiation of adherent NS cells. RHB-Basal does not contain any growth factors or neuronal supplements. Therefore, customers can tailor RHB-Basal to suit the specific requirements of their cell type by the addition of preferred supplements.

Storage: Upon receipt, store at -20°C until ready to use. Under these storage conditions, the product is stable for 12 months from the date of manufacture (see label).

Once thawed, store at 4°C and use within 4 weeks.

This product is light sensitive, and should be protected from light.

Preparation: Thaw the medium in a water bath (37°C) in the dark, and remove the medium from the water bath just before it has completely thawed (i.e., do not allow the medium to warm up). Then, mix gently and thaw completely. Alternatively, thaw the medium at 4°C while protecting from light. If a precipitate appears, leave the medium at 4°C overnight to completely dissolve the precipitate. Do not use medium with visible precipitate; ensure it is dissolved before use.

RHB-Basal is a base medium to which customers can add supplements of their choice to meet specific cell type requirements.

Quality Control: SC Proven™ media products undergo rigorous quality control procedures before release.

Note

This product is for research use only. It is not intended for use in therapeutic or diagnostic procedures for humans or animals. Also, do not use this product as food, cosmetic, or household item, etc. Takara products may not be resold or transferred, modified for resale or transfer, or used to manufacture commercial products without written approval from TAKARA BIO INC.

If you require licenses for other use, please contact us by phone at +81 77 565 6973 or from our website at www.takara-bio.com.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product web page. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

All trademarks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions.