



5-Fluoroorotic Acid

Cat. No. 10836-013

Size: 1 g

Store at -20°C.

Description: 5-Fluoroorotic acid (5FOA) inhibits cell growth in numerous organisms that utilize orotic acid as a source of the pyrimidine ring (1-8). 5FOA is converted to fluoroorotidine monophosphate and subsequently decarboxylated to form 5-fluorouridine monophosphate. Growth inhibition is presumably due to the formation of fluorodeoxyuridine, which is a potent inhibitor of thymidylate synthetase. Mutations in either of the two enzymes mediating these steps can be selected with 5FOA (9). In *Saccharomyces cerevisiae*, both *ura3⁻* and *ura5⁻* mutants can grow on plates containing 5FOA (1) whereas growth of cells expressing wild-type URA3 or URA5 is inhibited. In addition, events that decrease expression levels of URA3 can also be identified using 5FOA (10).

This material is recommended for use with the ProQuest™ Two-Hybrid System with Gateway™ Technology (Cat. No. 10835-031).

Caution: This product is considered hazardous. Please consult the Material Safety Data Sheet for health and safety information.

Quality Control: This product has been qualified in an inhibition assay using yeast control strains from the ProQuest™ Two-Hybrid System.

Doc. Rev. 082901

This product is distributed for laboratory research only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Invitrogen Tech-LineSM U.S.A. 800 955 6288

References:

1. Boeke, J.D., Trueheart, J., Natsoulis, G., and Fink G.R. (1987) *Methods Enzymol* 154, 164.
2. Krooth, R.S., Hsiao, W.L., and Potvin, B.W. (1979) *Somatic Cell Genet* 5, 551.
3. Worsham, P.L., and Goldman, W.E. (1988) *Mol Gen Genet* 214, 348.
4. Rathod, P.K., Khatri, A., Hubbert, T., and Milhous, W.K. (1989) *Antimicrob Agents Chemother* 33, 1090.
5. Gruber, F., Visser, J., Kubicek, C.P., and de Graaff, L.H. (1990) *Curr Genet* 18, 71.
6. Skory, C.D., Horng, J.S., Pestka, J.J., and Linz, J.E. (1990) *Appl Environ Microbiol* 56, 3315.
7. Kwon-Chung, K.J., Varma, A., Edman, J.C., and Bennett, J.E. (1992) *J Med Vet Mycol* 30, 61.
8. Yang, V.W., Marks, J.A., Davis, B.P., and Jeffries, T.W. (1994) *Appl Environ Microbiol* 60, 4245.
9. Kaiser, C., Michaelis, S., and Mitchell, A. (1994) *Methods in Yeast Genetics*. Cold Spring Harbor Laboratory Press.
10. Vidal, M., Brachmann, R.K., Fattaey, A., Harlow, E., and Boeke, J.D. (1996) *Proc. Natl. Acad. Sci. USA* 93, 10315.