

Yeast reference strains

Cat. #YSC1058, YSC1059, YSC1060, YSC1071, YSC1764

Storage

- 4 °C for up to one week
- -80 °C indefinitely

Product description

Culture of *Saccharomyces cerevisiae* in YPD broth with an inert growth indicator + 15% glycerol.

W303 MATa	MATa ura3-52 trp1Δ2 leu2-3_112 his3-11 ade2-1 can1-100
W303 MATα	MATα ura3-52 trp1Δ2 leu2-3_112 his3-11 ade2-1 can1-100
W303 Het-erozygous Diploid	MATa/MATα ura3-52/ura3-52 trp1Δ2/ trp1Δ2 leu2-3_112/ leu2-3_112 his3-11/ his3-11 ade2-1 can1-100/ ade2-1 can1-100
S288C	MATa p ^o trp1-0
Y800 ²	[MATa leu2-Δ98cry1 ^R /MATα leu2-Δ98CRY1 ade2-101 HIS3/ ade2-101 his3-Δ200 ura3-52 can1 ^R /ura3-52CAN1 lys2-801/ lys2-801 CYH2/cyh2 ^R trp1-1/TRP1 Cir ^R] carrying pGAL-cre (amp,ori, CEN, LEU2)

Useful websites and references

Saccharomyces Genome Deletion Project, click [here](#).

1. Link, J.L., and Olson M.V. Physical Map of the *Saccharomyces cerevisiae* Genome at 110-Kilobase Resolution. *Genetics* **127**, 681-698, 1991.
2. Burns, N., Grimwade, B., Ross-Macdonald, P.B., Choi, E.-Y., Finberg, K., Roeder, G.S., and Snyder, M. 1994. Large-scale characterization of gene expression, protein localization and gene disruption in *Saccharomyces cerevisiae*. *Genes & Dev.* **8**, 087-1105.

If you have any questions, contact

t +44 (0) 1223 976 000 (UK) **or** +1 800 235 9880 (USA); +1 303 604 9499 (USA)

f + 44 (0)1223 655 581

w horizondiscovery.com/contact-us **or** dharmacon.horizondiscovery.com/service-and-support

Horizon Discovery, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom

©2018 Horizon Discovery Group Company—All rights reserved. All trademarks are the property of Horizon Discovery Company unless otherwise specified. First published June 2014. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.