

# Yeast Tet-promoters Hughes (yTHC) parental strain—R1158

Cat. #YSC1210

The Yeast Tet-promoters Hughes (yTHC) Parental Strain contains 632 essential yeast genes for which expression is regulated by doxycycline. The endogenous promoter has been replaced with a TET-titratable promoter in the genome. Thus, the expression of the gene can be switched off by the addition of doxycycline to the yeast's growth medium.

The yTHC mutant strains are provided in the haploid MATa strain R1158. This strain was created from the background strain BY4741 by a one-step integration of the tTA transactivator, under the control of the CMV promoter, at the URA3 locus (*Hughes et al., Cell.* 2000 Jul 7;102(1):109-26).

## Storage

4 °C for up to one week

-80 °C indefinitely

## Product description

Culture of *S. cerevisiae* in YPD broth + 15% glycerol.

## Making a stock culture

Once the strain has been streaked, we recommend making a stock of the pure culture. Inoculate the pure culture in YPD broth and incubate for 48 hours at 30 °C. Transfer 850 µL of culture into a sterile polypropylene tube and add 150 µL sterile glycerol to make a 15% glycerol freezing solution. Vortex the culture to evenly mix the glycerol throughout the culture. The culture can be stored indefinitely at -80 °C.

## Genotype of *S. cerevisiae* R1158

URA::CMV-tTA MATa his3-1 leu2-0met15-0

## 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](http://horizondiscovery.com/contact-us) **or** [dharmacon.horizondiscovery.com/service-and-support](http://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 2015. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.