

Yeast Tet-promoters Hughes (yTHC) collection

Applicable catalog numbers

Cat. #	Description
YSC1182	Yeast Tet-promoters Hughes (yTHC) Collection
YSC1180	Yeast Tet-promoters Hughes (yTHC) Strain
YSC1210	Yeast Tet-promoters Hughes Parental Strain - R1158

Product description

The creation of a genome-wide [yeast deletion collection](#), reveals that there are over 1000 protein-coding genes that are essential for viability. The function of many of these essential genes has yet to be defined. In part this is due to the difficulty of utilizing essential genes in functional assays. The Tet-promoters Yeast Collection from the Hughes Laboratory, University of Toronto, helps to overcome this barrier.

The Yeast Tet-promoters Hughes (yTHC) Collection contains over 800 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.

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 2015. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.

Quality control

All cultures are checked for growth prior to shipment.

Shipping and storage

yTHC constructs are distributed as cultures of *S. cerevisiae* in YPD broth + G418 (300 g/mL) + 15% glycerol. Individual constructs are shipped at room temperature. Collections are shipped in 96-well plate format on dry ice. Individual constructs and collections should be stored at -80°C .

To allow any CO₂ that may have dissolved into the medium from the dry ice in shipping to dissipate, please store constructs at -80°C for at least 48 hours before thawing.