

CITATIONS

DUBprofiler – Oct 2022

The application of DUBprofiler™ service to evaluate the selectivity of USP30 inhibitors by Eisai

1: High-content phenotypic screen to identify small molecule enhancers of Parkin-dependent ubiquitination and mitophagy

Roberta Tufi, Emily H. Clark, Tamaki Hoshikawa, Christiana Tsagkaraki, Jack Stanley, Kunitoshi Takeda, James M. Staddon, Thomas Briston

bioRxiv preprint doi:
<https://doi.org/10.1101/2022.09.30.509127>

The application of DUBprofiler service in SARS-CoV-2 PLpro small molecule drug discovery by Walter and Eliza Hall Institute/Monash University/CSIRO/The University of Melbourne

2: Insights Into Drug Repurposing, as Well as Specificity and Compound Properties of Piperidine-Based SARS-CoV-2 PLpro Inhibitors

Dale J. Calleja, Nathan Kuchel, Bernadine G. C. Lu, Richard W. Birkinshaw, Theresa Klemm,

Marcel Doerflinger, James P. Cooney, Liana Mackiewicz, Amanda E. Au, Yu Q. Yap, Timothy R Blackmore, Kasiram Katneni, Elly Crighton, Janet Newman, Kate E. Jarman, Melissa J. Call, Bernhard C. Lechtenberg, Peter E. Czabotar, Marc Pellegrini, Susan A. Charman, Kym N. Lowes, Jeffrey P. Mitchell, Ueli Nachbur, Guillaume Lessene, David Komander

Front Chem. 2022; 10: 861209. Published online 2022 Apr 12. doi: 10.3389/fchem.2022.861209

PMCID: PMC9039177

The application of DUBprofiler service in identifying novel DUB substrates by the Dana-Farber Cancer Institute and Harvard Medical School

3: Proteomics-Based Identification of DUB Substrates Using Selective Inhibitors

Cell Chem Biol. Author manuscript; available in PMC 2022 Jan 21. Published in final edited form as: Cell Chem Biol. 2021 Jan 21; 28(1): 78–87.e3. Published online 2020 Oct 1. doi: 10.1016/j.chembiol.2020.09.005

PMCID: PMC7855594

The application of DUBprofiler service in identifying novel DUB substrates by University College London and Eisai

4: Investigation of USP30 inhibition to enhance Parkin-mediated mitophagy: tools and approaches

Eliona Tsefou, Alison S. Walker, Emily H. Clark, Amy R. Hicks, Christin Luft, Kunitoshi Takeda, Toru Watanabe, Bianca Ramazio, James M. Staddon, Thomas Briston, Robin Ketteler

Biochem J. 2021 Dec 10; 478(23): 4099–4118. Prepublished online 2021 Oct 27. Published online 2021 Dec 6. doi: 10.1042/BCJ20210508

The application of DUBprofiler service in identifying novel DUB substrates by the University of Liverpool and University of Sheffield

5: Benchmarking a highly selective USP30 inhibitor for enhancement of mitophagy and pexophagy

Emma V Rusilowicz-Jones, Francesco G Barone, Fernanda Martins Lopes, Elezabeth Stephen, Heather Mortiboys, Sylvie Urbé, Michael J Clague

Life Sci Alliance. 2022 Feb; 5(2): e202101287. Published online 2021 Nov 29. doi: 10.26508/lsa.202101287

PMCID: PMC8645336

The application of DUBprofiler service in identifying novel DUB substrates by the University of Liverpool, University of Oxford, FORMA Therapeutics, Max-Planck-Institute of Molecular Physiology, Technische Universität

Dortmund, University of Sheffield, Walter and Eliza Hall Institute of Medical Research, University of Melbourne and University of Newcastle

6: USP30 sets a trigger threshold for PINK1–PARKIN amplification of mitochondrial ubiquitylation

Emma V Rusilowicz-Jones, Jane Jardine, Andreas Kallinos, Adan Pinto-Fernandez, Franziska Guenther, Mariacarmela Giurrandino, Francesco G Barone, Katy McCarron, Christopher J Burke, Alejandro Murad, Aitor Martinez, Elena Marcassa, Malte Gersch, Alexandre J Buckmelter, Katherine J Kayser-Bricker, Frederic Lamoliatte, Akshada Gajbhiye, Simon Davis, Hannah C Scott, Emma Murphy, Katherine England, Heather Mortiboys, David Komander, Matthias Trost, Benedikt M Kessler, Stephanos Ioannidis, Michael K Ahljanian, Sylvie Urbé, Michael J Clague

Life Sci Alliance. 2020 Aug; 3(8): e202000768. Published online 2020 Jul 7. doi: 10.26508/lsa.202000768

PMCID: PMC7362391

The application of DUBprofiler service in DUB targeting small molecule drug discovery by the Dana-Farber Cancer Institute and Harvard Medical School

7: Selective USP7 inhibition elicits cancer cell killing through a p53-dependent mechanism

Nathan J. Schauer, Xiaoxi Liu, Robert S. Magin, Laura M. Doherty, Wai Cheung Chan, Scott B. Ficarro, Wanyi Hu, Rebekka M. Roberts, Roxana E. Jacob, Björn Stolte, Andrew O. Giacomelli, Sumner Perera, Kyle McKay, Sarah A. Boswell, Ellen L. Weisberg, Arghya Ray, Dharminder Chauhan, Sirano Dhe-Paganon, Ken C. Anderson, James D. Griffin, Jianing Li, William C.

Hahn, Peter K. Sorger, John R. Engen, Kimberly Stegmaier, Jarrod A. Marto, Sara J. Buhrlage

Sci Rep. 2020; 10: 5324. Published online 2020 Mar 24. doi: 10.1038/s41598-020-62076-x
PMCID: PMC7093416

The application of DUBprofiler service in DUB targeting small molecule drug discovery by Almac Discovery and Queen's University

8: Identification and Structure-Guided Development of Pyrimidinone Based USP7 Inhibitors

Colin R. O'Dowd, Matthew D. Helm, J. S. Shane Rountree, Jakub T. Flasz, Elias Arkoudis, Hugues Miel, Peter R. Hewitt, Linda Jordan, Oliver Barker, Caroline Hughes, Ewelina Rozycka, Eamon Cassidy, Keeva McClelland, Ewa Odrzywol, Natalie Page, Stephanie Feutren-Burton, Scarlett Dvorkin, Gerald Gavory, Timothy Harrison

ACS Med Chem Lett. 2018 Mar 8; 9(3): 238–243. Published online 2018 Feb 21. doi: 10.1021/acsmchemlett.7b00512

PMCID: PMC5846043

The application of DUBprofiler service in identifying novel DUB substrates by the Mitobridge, Inc. (an Astellas company) and Aurigene Discovery Technologies

9: Novel highly selective inhibitors of ubiquitin specific protease 30 (USP30) accelerate mitophagy

Arthur F Kluge 1, Bharat R Lagu 2, Pranab Maiti 3, Mahaboobi Jaleel 3, Michael Webb 1, Jyoti Malhotra 1, Ashley Mallat 1, P Akhila Srinivas 3, James E Thompson 1

Bioorg Med Chem Lett 2018 Aug 15;28(15):2655–2659. doi: 10.1016/j.bmcl.2018.05.013. Epub 2018 May 8

The application of DUBprofiler service in DUB targeting small molecule drug discovery by Almac Discovery and Queen's University

10: Discovery and characterization of highly potent and selective allosteric USP7 inhibitors.

Gerald Gavory, Colin R O'Dowd, Matthew D Helm, Jakub Flasz, Elias Arkoudis, Anthony Dossang, Caroline Hughes, Eamon Cassidy, Keeva McClelland, Ewa Odrzywol, Natalie Page, Oliver Barker, Hugues Miel & Timothy Harrison

Nature Chemical Biology volume 14, pages118–125 (2018) and (2017) Nat Chem Biol. doi: 10.1038/nchembio.2528. [Epub ahead of print]

The application of DUBprofiler service in DUB targeting small molecule drug discovery by the Dana-Farber Cancer Institute and Harvard Medical School

11: Structure-guided development of a potent and selective noncovalent active site inhibitor of USP7

Ilaria Lamberto, Xiaoxi Liu, Hyuk-Soo Seo, Nathan J Schauer, Roxana E Iacob, Wanyi Hu, Deepika Das, Tatiana Mikhailova, Ellen L Weisberg, John R Engen, Kenneth C Anderson, Dharminder Chauhan, Sirano Dhe-Paganon, Sara J Buhrlage

Cell Chem Biol. Author manuscript; available in PMC 2018 Dec 21. Published in final edited form as: Cell Chem Biol. 2017 Dec 21; 24(12): 1490–1500.e11. Published online 2017 Oct 19. doi: 10.1016/j.chembiol.2017.09.003

PMCID: PMC5749250

The application of DUBprofiler service in DUB targeting small molecule drug discovery by the University of Pittsburgh

12: Targeting the deubiquitinase STAMBP inhibits NALP7 inflammasome activity

Joseph S. Bednash, Nathaniel Weathington, James Londino, Mauricio Rojas, Dexter L. Gulick, Robert Fort, SeungHye Han, Alison C. McKelvey, Bill B. Chen, Rama K. Mallampalli

Nat Commun. 2017; 8: 15203. Published online 2017 May 11. doi: 10.1038/ncomms15203

PMCID: PMC5437278

The application of DUBprofiler service in DUB targeting small molecule drug discovery by Genentech

13: USP7 small-molecule inhibitors interfere with ubiquitin binding.

Lorna Kategaya 1 2, Paola Di Lello 3, Lionel Rougé 3, Richard Pastor 4, Kevin R Clark 5, Jason Drummond 5, Tracy Kleinheinz 5, Eva Lin 1, John-Paul Upton 1 2, Sumit Prakash 1 2, Johanna Heideker 1 2, Mark McClelland 1 2, Maria Stella Ritorto 6, Dario R Alessi 6, Matthias Trost 7, Travis W Bainbridge 8, Michael C M Kwok 8, Taylur P Ma 9, Zachary Stiffler 10, Bradley Brasher 10, Yinyan Tang 11, Priyadarshini Jaishankar 11, Brian R Hearn 11, Adam R Renslo 11, Michelle R Arkin 11, Frederick Cohen 4, Kebin Yu 9, Frank Peale 12, Florian Gnad 13, Matthew T Chang 13, Christiaan Klijn 13, Elizabeth Blackwood 14, Scott E Martin 1, William F Forrest 13, James A Ernst 8, Chudi Ndubaku 4, Xiaojing Wang 4, Maureen H Beresini 5, Vickie Tsui 4, Carsten Schwerdtfeger 10, Robert A Blake 5, Jeremy Murray 3, Till Maurer 3, Ingrid E Wertz Kategaya et al. (2017)

Nature. 550(7677):534-538. doi: 10.1038/nature24006. Epub 2017 Oct 18.

The application of DUBprofiler service in DUB targeting small molecule drug discovery by Progenra, University of Connecticut and Yale University

14: USP7-Specific Inhibitors Target and Modify the Enzyme's Active Site via Distinct Chemical Mechanisms.

Alexandra Pozhidaeva 1, Gabrielle Valles 1, Feng Wang 2, Jian Wu 2, David E Sterner 2, Phuong Nguyen 2, Joseph Weinstock 2, K G Suresh Kumar 2, Jean Kanyo 3, Dennis Wright 4, Irina Bezsonova 5

Cell Chem Biol. 2017 Dec 21;24(12):1501-1512.e5. doi: 10.1016/j.chembiol.2017.09.004. Epub 2017 Oct 19.

The application of DUBprofiler service in DUB targeting small molecule drug discovery by FORMA Therapeutics

15: Molecular basis of USP7 inhibition by selective small-molecule inhibitors

Andrew P. Turnbull, Stephanos Ioannidis, Wojciech W. Krajewski, Adan Pinto-Fernandez, Claire Heride, Agnes C.L. Martin, Louise M. Tonkin, Elizabeth C. Townsend, Shane M. Buker, David R. Lancia, Jr, Justin A. Caravella, Angela V. Toms, Thomas M. Charlton, Johanna Lahdenranta, Erik Wilker, Bruce C. Follows, Nicola J. Evans, Lucy Stead, Cristina Alli, Vladislav V. Zarayskiy, Adam C. Talbot, Alexandre J. Buckmelter, Minghua Wang, Crystal L. McKinnon, Fabienne Saab, Joanna F. McGouran, Hannah Century, Malte Gersch, Marc S. Pittman, C. Gary Marshall, Tony M. Raynham, Mary Simcox, Lorna M.D. Stewart, Sheila B. McLoughlin, Jaime A. Escobedo, Kenneth W.

Bair, Christopher J. Dinsmore, Tim R. Hammonds, Sunkyu Kim, Sylvie Urbé, Michael J. Clague, Benedikt M. Kessler, David Komander

Nature. Author manuscript; available in PMC 2018 Jul 3. Published in final edited form as: Nature. 2017 Oct 26; 550(7677): 481–486. Published online 2017 Oct 18. doi: 10.1038/nature24451

PMCID: PMC6029662

The application of DUBprofiler service in DUB targeting small molecule drug discovery by Almac Discovery and Astra Zeneca

16: Identification and characterisation of dual inhibitors of the USP25/28 Deubiquitinating enzyme subfamily.

Jonathan D Wrigley 1, Gerald Gavory 2, Iain Simpson 3, Marian Preston 1, Helen Plant 1, Jenna Bradley 1, Anne U Goeppert 1, Ewelina Rozycka 2, Gareth Davies 1, Jarrod Walsh 1, Andrea Valentine 2, Keeva McClelland 2, Krzysztofa Ewa Odrzywol 2, Jonathan Renshaw 1, Joanna Boros 1, Jonathan Tart 1, Lindsey Leach 1, Thorsten Nowak 3, Richard A Ward 3, Timothy Harrison 2, David M Andrews 3 Wrigley et al. (2017)

ACS Chem Biol DOI: 10.1021/acscchembio.7b00334. Epub 2017 Nov 28.

The application of DUBprofiler service in identifying novel DUB substrates by the Linköping University, Mayo Clinic, Karolinska Institute and Uppsala University

17: Synthesis and Evaluation of Derivatives of the Proteasome Deubiquitinase Inhibitor b-AP15

Xin Wang, Pádraig D'Arcy, Thomas R. Caulfield, Aneel Paulus, Kasyapa Chitta, Chitralekha Mohanty, Joachim Gullbo, Asher Chanan-Khan, Stig Linder

Chem Biol Drug Des. Author manuscript; available in PMC 2016 Nov 1. Published in final edited form as: Chem Biol Drug Des. 2015 Nov; 86(5): 1036–1048. Published online 2015 May 27. doi: 10.1111/cbdd.12571

The application of Ubiquigent's DUBprofiler service in identifying novel DUB substrates by the Heinrich-Heine-University, The Norwegian Radium Hospital and University of Oslo

18: Deubiquitinase inhibition by WP1130 leads to ULK1 aggregation and blockade of autophagy

Stefan Drießen, Niklas Berleth, Olena Friesen, Antje S Löffler, Philip Böhler, Nora Hieke, Fabian Stuhldreier, Christoph Peter, Kay O Schink, Sebastian W Schultz, Harald Stenmark, Petter Holland, Anne Simonsen, Sebastian Wesselborg, Björn Stork

Autophagy. 2015 Sep; 11(9): 1458–1470. Published online 2015 Jul 24. doi: 10.1080/15548627.2015.1067359

PMCID: PMC4590650

Please let us know if you have used *DUBprofiler* in your publication so that we can keep our citation list up to date. Thank you!

Email: services@ubiquigent.com