

CITATIONS

UBISITE*profiler* – May 2023

1: UbiSite approach for comprehensive mapping of lysine and N-terminal ubiquitination sites.

Akimov V, Barrio-Hernandez I, Hansen SVF, Hallenborg P, Pedersen AK, Bekker-Jensen DB, Puglia M, Christensen SDK, Vanselow JT, Nielsen MM, Kratchmarova I, Kelstrup CD, Olsen JV, Blagoev B.

Nat Struct Mol Biol. 2018 Jul;25(7):631-640. doi: 10.1038/s41594-018-0084-y. Epub 2018 Jul 2. PMID: 29967540.

2: Deubiquitinating enzymes and the proteasome regulate preferential sets of ubiquitin substrates.

Trulsson F, Akimov V, Robu M, van Overbeek N, Berrocal DAP, Shah RG, Cox J, Shah GM, Blagoev B, Vertegaal ACO.

Nat Commun. 2022 May 18;13(1):2736. doi: 10.1038/s41467-022-30376-7. PMID: 35585066; PMCID: PMC9117253.

3: Identification of ester-linked ubiquitylation sites during TLR7 signalling increases the

number of inter-ubiquitin linkages from 8 to 12.

McCrary EH, Akimov V, Cohen P, Blagoev B.

Biochem J. 2022 Dec 9;479(23):2419-2431. doi: 10.1042/BCJ20220510. PMID: 36408944; PMCID: PMC9788571.

4: Large-Scale Profiling of Unexpected Tryptic Cleaved Sites at Ubiquitinated Lysines.

Sun Z, Xiao W, Li N, Chang L, Xu P, Li Y.

J Proteome Res. 2023 Apr 7;22(4):1245-1254. doi: 10.1021/acs.jproteome.2c00748. Epub 2023 Mar 6. PMID: 36877145.



Please let us know if you use **UBISITEprofiler™** in your research so that we can keep our citation list up to date. Thank you!

Email: services@ubiquigent.com