

Simultaneous targeted and discovery phosphoproteomics of cell signaling pathways using novel hybrid-DIA acquisition strategy

Ana Martínez-Val¹, Kyle Fort², Giulia Franciosa¹, Claire Koenig¹, Daniel Lopez Ferrer³, Andreas F. Huhmer³, Yasushi Ishihama⁴, Yu-ju Chen⁵, Yue Xuan², Alexander Makarov², Jesper V. Olsen¹

¹Novo Nordisk Foundation Center for Protein Research, University of Copenhagen, DENMARK

³Thermo Fisher Scientific, San Jose, USA

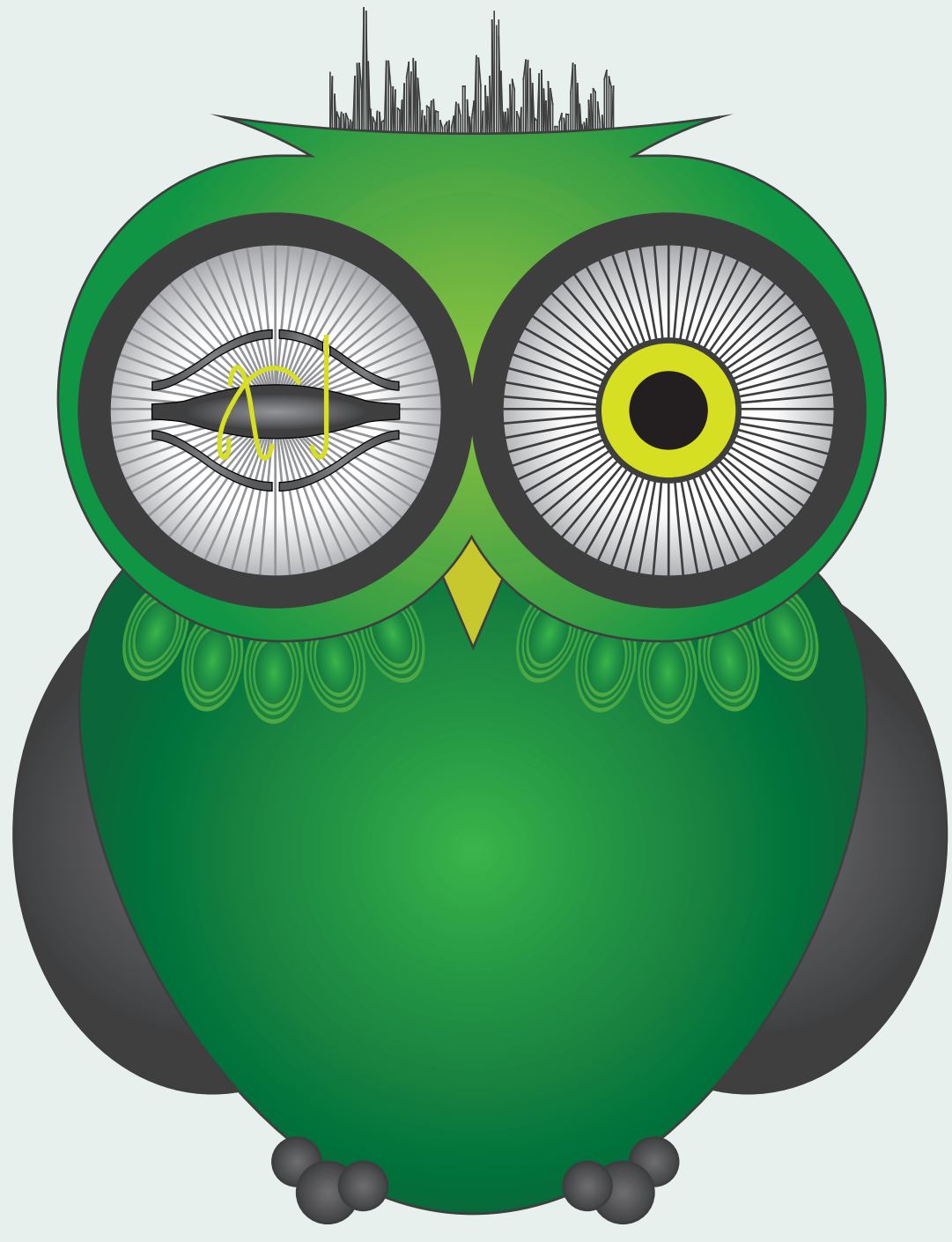
⁵Academia Sinica, Taipei, TAIWAN

²Thermo Fisher Scientific, Bremen, GERMANY

⁴Kyoto University, Kyoto, JAPAN

ana.mdval@cpr.ku.dk

OLSEN
GROUP

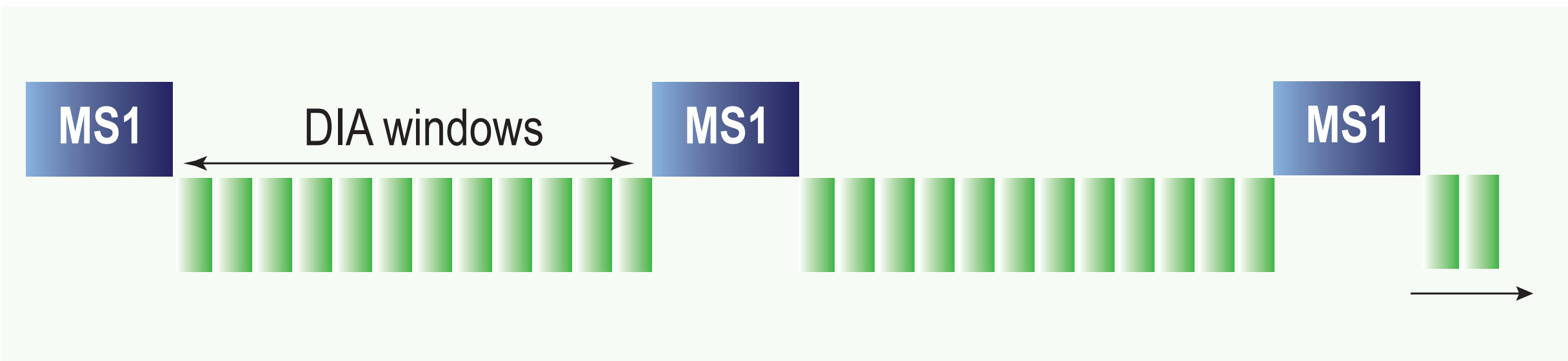


Aims

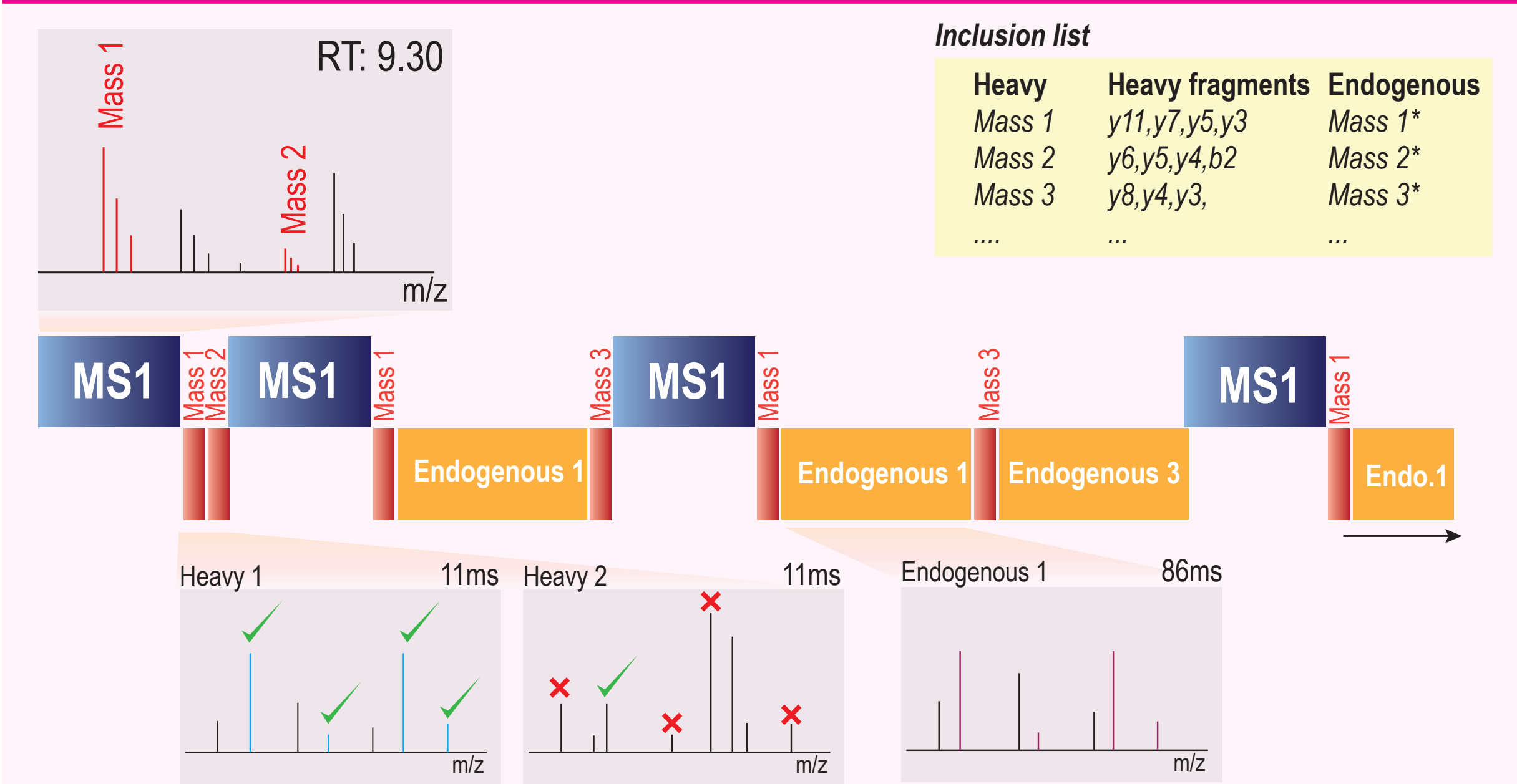
- **Hybrid DIA** is based on an **Application Programming Interface**, which enables the ability to **combine targeted and discovery proteomics**, by intercalating DIA scans with accurate triggering of predefined peptide targets.
- We have **benchmarked the sensitivity** of hybrid DIA against purely targeted methods, such as SureQuant.
- We have evaluated the utility of this data acquisition strategy to study changes in phosphorylation levels of 131 relevant peptides together with a comprehensive analysis of the phospho-proteome.

Overview of Data Acquisition Methods

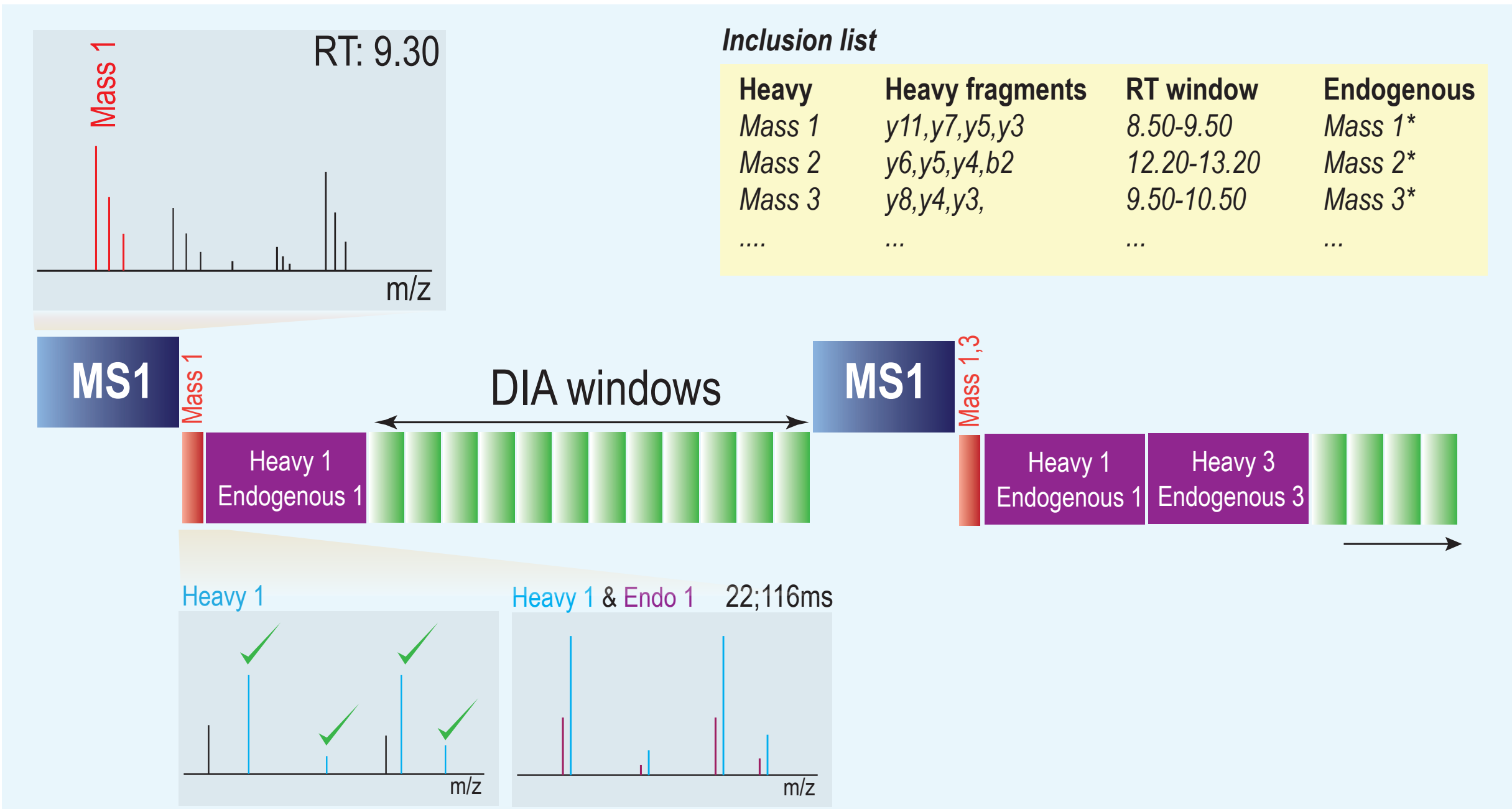
Conventional DIA



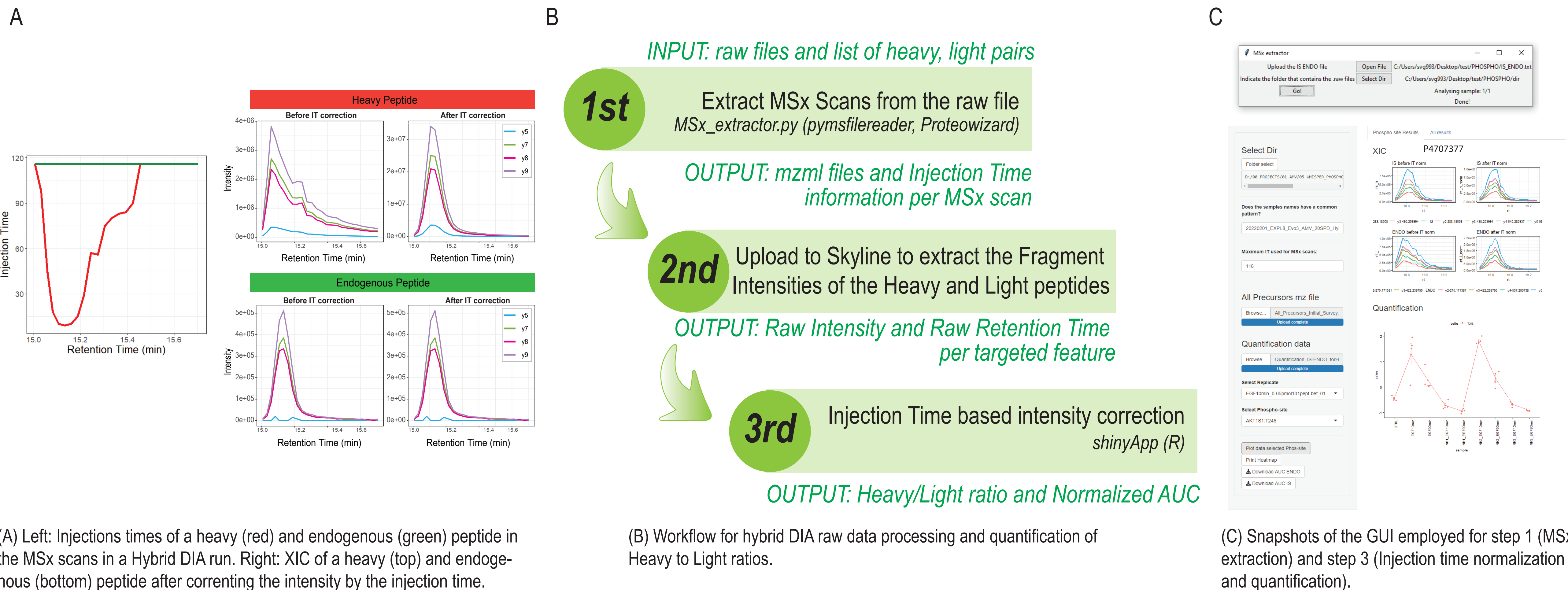
Targeted method (i.e. SureQuant)



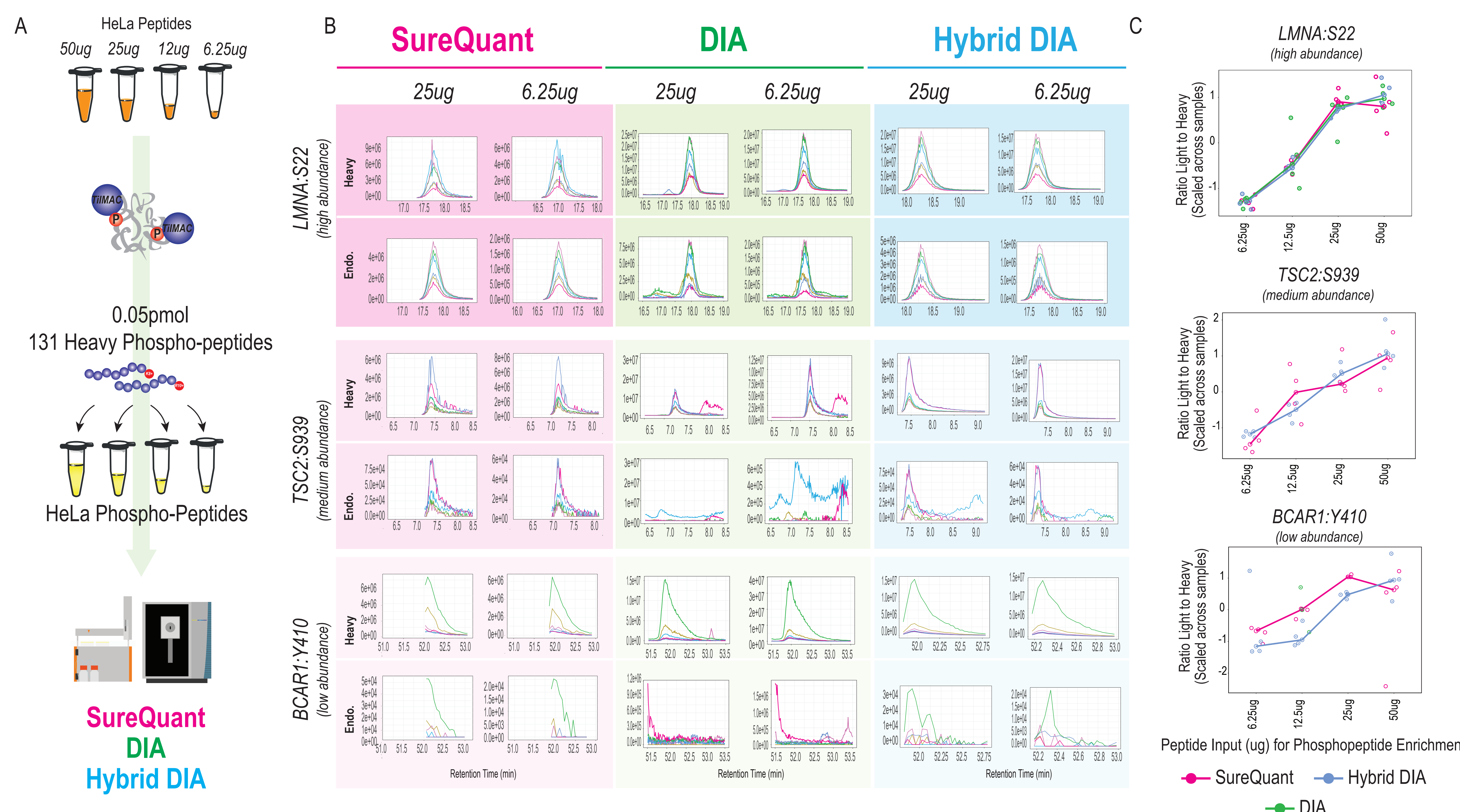
Hybrid DIA



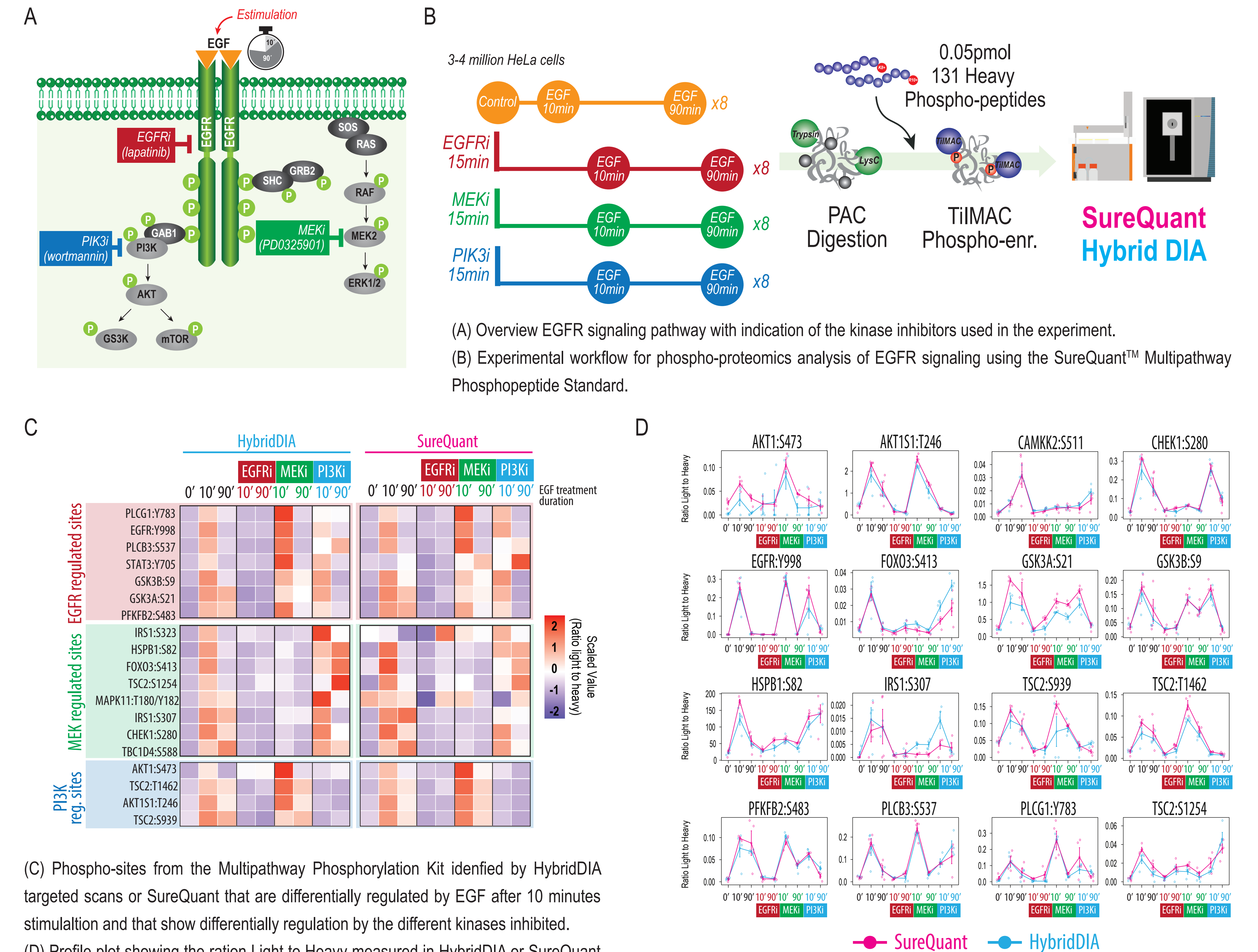
Data Analysis Pipeline



Improved limit of detection



Targeted and discovery analysis of EGF stimulation



(C) Phospho-sites from the Multipathway Phosphorylation Kit identified by HybridDIA targeted scans or SureQuant that are differentially regulated by EGF after 10 minutes stimulation and that show differentially regulation by the different kinases inhibited.

(D) Profile plot showing the ration Light to Heavy measured in HybridDIA or SureQuant for some relevant EGFR targets.

E

**Full phosphoproteome analysis:
Spectronaut directDIA results**

5,055 Phospho-sites identified (Loc. Prob >0.75)

3,611 Phospho-sites identified quantified in at least 75% of the replicates.

1,214 Phospho-sites differential in total (ANOVA FDR<0.05%)

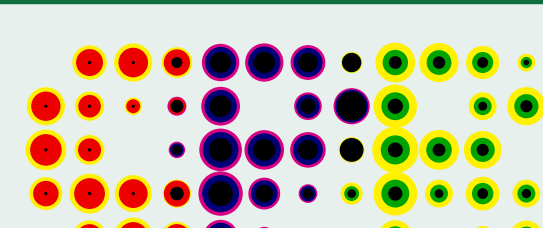
224 Phospho-sites activated due to EGF stimulation (10 minutes)

(E) Full phosphoproteome results obtained from a library-free search of the HybridDIA files.

(F) Kinase activity inferred from the full phospho-proteome profiles in the different experimental condition calculated with RoKAI App (v2.1.3).

Acknowledgements: **ThermoFisher**
SCIENTIFIC

Funding: NNF14CC0001
EPIC-XS-823839



Novo Nordisk Foundation
Center for Protein Research