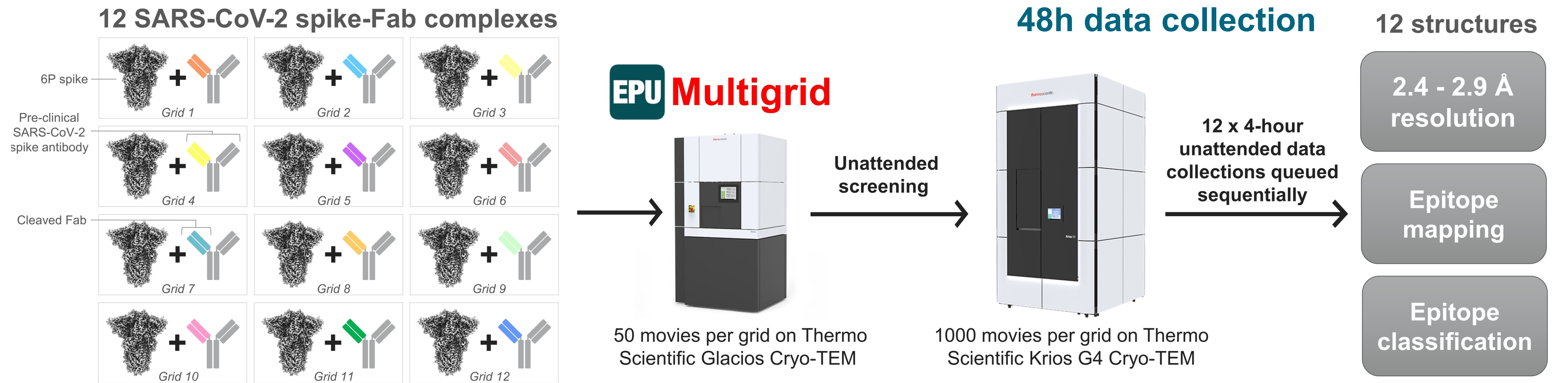


# Cryo-EM: High speed epitope mapping

Ieva Drulyte<sup>1</sup>, Stefan Koester<sup>2</sup>, DiannaLundberg<sup>2</sup>, Kurt Eng<sup>2</sup>, Andrey Malyutin<sup>3</sup>, Aaron McGrath<sup>3</sup>, Chunyan Wang<sup>4</sup>, Frank van Kuppeveld<sup>4</sup>, Berend-Jan Bosch<sup>4</sup>, Mazdak Radjainia<sup>1</sup>, Jerry Thomas<sup>3</sup>, Maureen Magnay<sup>2</sup> and Daniel Hurdiss<sup>4</sup>

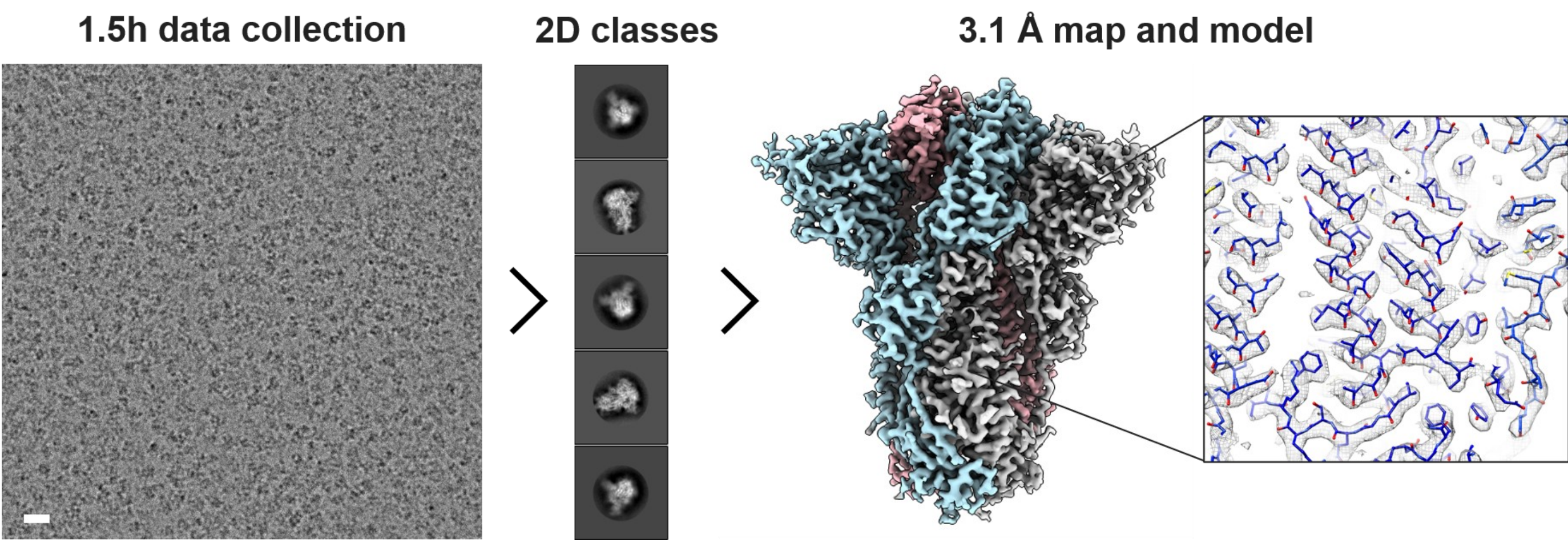
<sup>1</sup>Materials and Structural Analysis Division, Thermo Fisher Scientific, Netherlands; <sup>2</sup>Takeda Pharmaceuticals, Cambridge MA, USA; <sup>3</sup> Takeda Pharmaceuticals, San Diego, CA, USA; <sup>4</sup> Department of Biomolecular Health Sciences, Utrecht University, Netherlands.

## 12 structures in 48h



Cryo-electron microscopy (cryo-EM) is a powerful tool for epitope mapping of antibodies that block SARS-CoV-2 virus entry. Defining the epitopes of neutralizing antibodies allows us to understand how antibodies can confer protective immunity against SARS-CoV-2. Here we show how 12 sub-3-angstrom reconstructions of spike protein in complex with 12 distinct Fabs can be obtained from a 2-day unattended microscopy session using EPU Multigrid.

METHODS



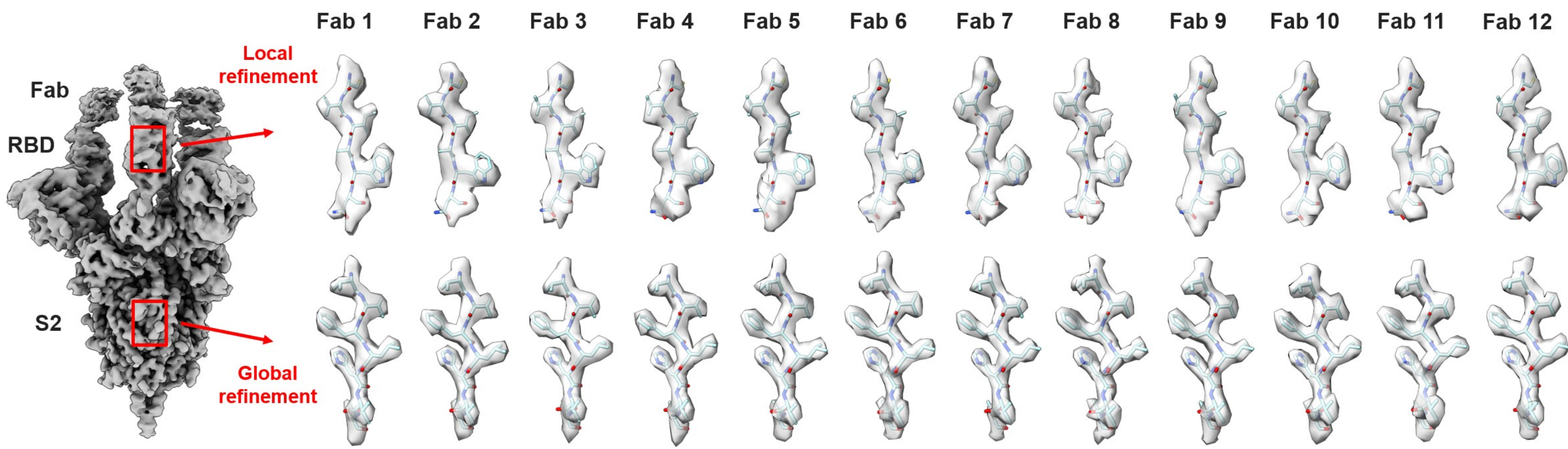
| Grid                   | Quantifoil R1.2/1.3                             |
|------------------------|---|
| Camera                 | Selectris X Energy Filter and Falcon 4 Detector |
| Slit width (eV)        | 10  |
| Nominal magnification  | 130,000x  |
| Pixel size (Å)         | 0.929   |
| Dose rate (e-/pix/sec) | 5.3   |
| Exposure time (sec)    | 8.32  |
| Total dose (e-/Å²)     | 48  |
| Fractionation          | EER   |
| Autofocus              | After centering                                 |
| Hole centering         | AFIS (6 µm image shift)                         |
| Number of images       | 12 x 1000                                       |
| Throughput             | ~250/hour                                       |

Table 1. Cryo-EM data acquisition parameters.

## Methods

Spike-Fab-complex grids were vitrified in duplicate and screened using Glacios cryo-TEM. Twelve best grids, one grid per complex, were taken forward for a 48-hour unattended EPU Multigrid session using Krios G4 cryo-TEM.

RESULTS



| Spike-Fab complex | Global resolution (Å) | Local resolution (Å) |
|-------------------|-----------------------|----------------------|
| Fab 1             | 2.8                   | 3.7                  |
| Fab 2             | 2.7                   | 4.0                  |
| Fab 3             | 2.9                   | 3.7                  |
| Fab 4             | 2.9                   | 3.7                  |
| Fab 5             | 2.8                   | 4.2                  |
| Fab 6             | 2.5                   | 3.7                  |
| Fab 7             | 2.9                   | 3.4                  |
| Fab 8             | 2.4                   | 3.1                  |
| Fab 9             | 2.8                   | 3.9                  |
| Fab 10            | 2.6                   | 3.4                  |
| Fab 11            | 2.5                   | 3.4                  |
| Fab 12            | 2.9                   | 3.4                  |

Table 2. Global and local resolutions achieved for each spike-FAB complex

## Results

We have obtained 12 sub-3Å structures of the SARS-CoV-2 spike protein in complex with RBD-binding Fabs. Our reconstructions are well resolved at the epitope-paratope interface despite the relatively small size of each dataset. Sample optimization and screening, as well as the use of both the Selectris X Energy Filter and the E-CFEG resulted in high quality datasets.

## Conclusion

Significant advances have been made in the speed, quality, and automation of cryo-EM data collection. The cryo-EM's resolution revolution will be followed by a throughput revolution, which increases efficiency and reduces cost per dataset, and enable fragment-based drug discovery and epitope mapping of large antibody panels.

In summary, cryo-EM is now poised to emulate the success of macromolecular crystallography thanks to faster data collection and EPU Multigrid.

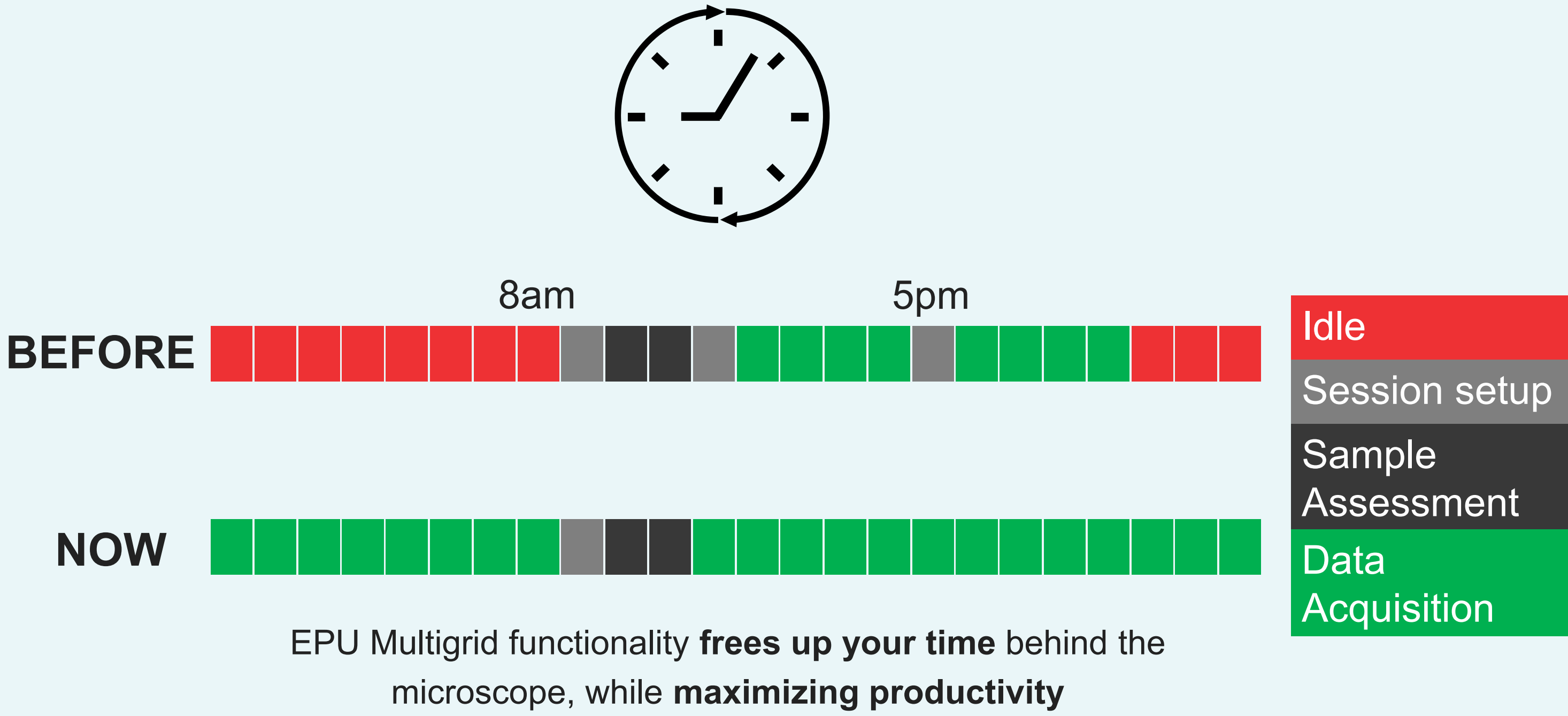
## Acknowledgements

This work was partially funded by the Corona Accelerated R&D in Europe (CARE) project. The CARE project has received funding from the Innovative Medicines Initiative 2 Joint Undertaking (JU) under grant agreement No 101005077. The JU receives support from the European Union's Horizon 2020 Research and Innovation Programme, the European Federation of Pharmaceutical Industries and Associations, the Bill & Melinda Gates Foundation, the Global Health Drug Discovery Institute, and the University of Dundee. The content of this publication only reflects the author's view and the JU is not responsible for any use that may be made of the information it contains. Antibody discovery was performed at Takeda Pharmaceuticals



ThermoFisher  
SCIENTIFIC

## Round-the-clock productivity and sample throughput



- 1

Acquire an Atlas on every grid to inspect samples
- 2

For each grid that you want to acquire:
  - Load and select Grid squares
  - Set foil hole settings such as ice filter and define an acquisition template. EPU will select the holes and prepare the grid squares automatically
  - Reload preferences or re-use previous session settings to set-up each grid in less than 15 minutes
- 3

Start a queued session

## Advantage of EPU Multigrid

- Reduce dedicated operator time from 8 to 3 hours for Glacios and Krios
- Fit sample assessment and high-resolution data acquisition in one session
- Tools can be used 24/7 due to unattended operation