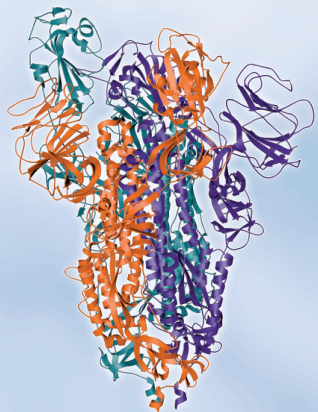


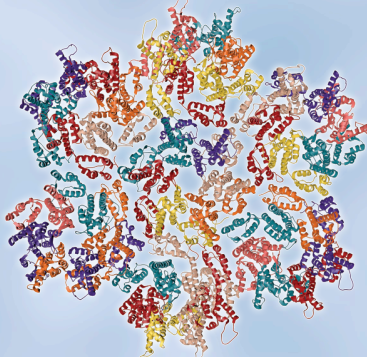
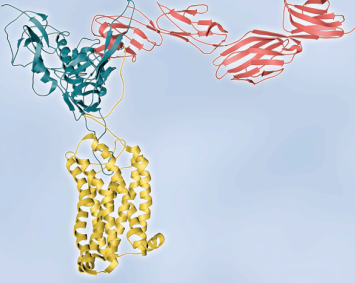



VIROLOGY		
SARS-COV-2		
First Spike Protein Structure, PDB 6VSB	Monoclonal Antibodies Binding Spike RBD, PDB 6XDG	3a Ion Channel, PDB 7KJR
		
SARS-CoV-2 binds to host cells via a trimeric spike glycoprotein. Researchers produced the first 3D structure of the multimeric protein by cryo-EM in mid-February 2020. The determination of the spike protein atomic structure was essential for developing diagnostics, antibody therapeutics, and vaccines. ²	Treatment with a single monoclonal antibody may create selective pressure for resistant mutants, so researchers are testing the efficacy of antibody cocktails that interact with the spike protein simultaneously. The cryo-EM structure showed REGN10933 (orange and purple) and REGN10987 (red and pink) antibodies binding the spike protein receptor-binding domain (RBD; turquoise). ³	Researchers have implicated the SARS 3a ion channel in processes such as viral release ⁴ and cell death. ⁵ In mice with a SARS 3a deletion mutation, viral titer and morbidity are reduced, ⁶ making 3a an effective vaccine and therapeutic target. Traditional structural analysis techniques are difficult to perform on membrane proteins such as ion channels; however, cryo-EM allowed researchers to view the SARS-CoV-2 3a dimer in its native conformation. ⁷

HUMAN IMMUNODEFICIENCY VIRUS (HIV-1)		
HIV-1 Capsid, PDB 3J34	HIV-1 gp120 Bound to Host Receptors, PDB 6MET	HIV-1 Envelope Glycan Shield, PDB 6X9R
		
X-ray crystallography determined only the structure of assembled HIV-1 capsid protein hexamers. Cryo-EM showed the structure and distribution of penton regions that facilitate capsid curvature and genome enclosing, and residues important for assembly and infectivity. ⁷	The HIV-1 envelope fusion protein gp120 (turquoise) interacts with host CD4 (pink) and a co-receptor such as CCR5 (yellow). Cryo-EM helped scientists understand the conformational changes that occur during these interactions and the role of CCR5 in fusing the virus to the host membrane. ⁹	The HIV-1 envelope (Env; yellow and red) is surrounded by a dense array of N-linked glycans (turquoise and green) that mask the protein from immune recognition. Using cryo-EM, researchers localized this “glycan shield” and discovered features of its structure and function that will aid in vaccine development. ¹⁰

SOLVING STRUCTURES IN A FLASH

SCIENTIFIC BREAKTHROUGHS WITH CRYOGENIC ELECTRON MICROSCOPY

Dramatic improvements to cryogenic electron microscopy (cryo-EM) technology have led to an explosion of research using this technique in diverse research areas. As cryo-EM protein structures increase in resolution, scientists learn more about molecular interactions than ever before, revolutionizing therapeutic development in virology, neurology, cancer, and a range of other diseases. Cryo-EM reveals key insights into the function, structure, and molecular basis for disease-associated mutations.




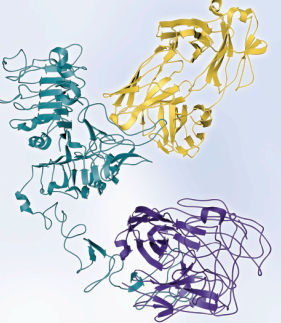
Traditional electron microscopy (TEM) lagged behind X-ray crystallography as the main choice for protein structure research. Although protein crystallization is challenging and time-consuming, traditional TEM techniques such as negative staining limit the determination of proteins to low-resolution structures that do not resolve the internal protein conformation, the interaction of amino acid side chains, and the atomic structure. That changed once Jacques Dubochet developed a method to vitrify (flash-freeze) electron microscopy samples; this approach prevents dehydration, avoids ice crystal formation, and captures proteins in their native state.¹ Vitrified cryo-EM specimens are illuminated with high-energy electrons (100-300 keV) and individual protein molecules are visualized in captured raw images. From these imaged molecules, scientists reconstruct high-resolution 3D maps and build atomic models of the protein structure.

Compared to X-ray crystallography, cryo-EM

- Does not require protein crystals
- Uses small amounts of purified protein
- Captures multiple protein conformations in their native state





NEUROSCIENCE	
C-terminal Portion of LRRK2, PDB 6VNO	Hallucinogen Complexed with a GPCR, PDB 6WHA
	
Leucine-rich repeat kinase 2 (LRRK2) protein is linked to genetically-inherited Parkinson's disease, but its mechanism and structure have been unknown for almost 20 years. Researchers finally visualized the protein in its natural environment with cryo-EM ¹¹ and cryo-electron tomography (cryo-ET), ¹² providing information that is crucial for future kinase inhibitor drug design.	Psychedelics are used as therapeutics for neuropsychiatric disorders such as depression, anxiety, and substance abuse. ¹³ These drugs activate the 5-HT _{2A} serotonin receptor (HTR2A)—a membrane G protein-coupled receptor (GPCR). To understand the mechanism of action, researchers performed cryo-EM to obtain the structure of a prototypical hallucinogen ligand (yellow) complexed within HTR2A (pink) and coupled G proteins (not shown). ¹⁴
CANCER	
Human CAK in Complex with Small Molecule Inhibitor, PDB 6XD3	Monoclonal Antibodies Simultaneously Binding to HER2, PDB 6GOE
	
Human CDK-activating kinase (CAK) is an important regulator of transcription initiation and the cell cycle. When CAK misregulates gene expression and cell division, cancer can occur, making it an attractive target for cancer treatment. ¹⁵ Using cryo-EM, scientists viewed how a small molecule inhibitor (yellow) binds to the human CAK complex (purple, turquoise, orange). ¹⁶	Human epidermal growth factor receptor 2 (HER2) associates with aggressive breast cancers and cancer recurrence. Interestingly, HER2-positive tumors are more likely to respond to treatment, especially combination antibody therapy. ¹⁷ To understand the synergy between two monoclonal antibodies, scientists used cryo-EM to image the simultaneous binding of trastuzumab (purple) and pertuzumab (yellow) antibodies to different HER2 subdomains (turquoise). ¹⁸

THE HISTORY OF CRYO-EM ^{19–22}	
1968	First 3D EM reconstruction; DeRosier and Klug
1974	First cryo-EM diffraction patterns; Taylor and Glaeser
1982	Vitrification developed; Dubochet et al.
1987	Birth of single particle reconstruction; van Heel, Radermacher, and Frank et al.
2007	Maximum-likelihood 3D software developed; Scheres et al.
2012	Direct electron-detection cameras first used; Grigorieff et al.
2015	Single particle cryo-EM named Method of the Year by <i>Nature Methods</i>
2017	Jacques Dubochet, Joachim Frank, and Richard Henderson win the Nobel Prize in Chemistry for cryo-EM
2020	Atomic level resolution achieved; Yip et al., Nakane et al.
2021 and beyond	Cryo-EM is poised to become the dominant technique for structure determination. New cryo-EM platforms are increasing ease-of-use and access.

Getting started in cryo-EM



including our single particle analysis (SPA) workflow, via step-by-step instructions and expert demonstrations.



The **Scientific Workflows App** (available in the iOS App Store and Google Play Store) puts the Cryo-EM University videos and content at your fingertips. The lite version (free) includes a lab notebook, lists of consumables and accessories, plus all of the rich useful content available at em-learning.com. The full version (part of the Accelerate service offering) also includes updated research articles every quarter, troubleshooting tips, and supervisor information.



Practical paths to accessing cryo-EM

Thermo Fisher Scientific provides accessible ways to leverage the advantages of cryo-EM without making a full investment. Our Cryo-EM Access program connects biopharma to world-class facilities where you can advance your research by obtaining cryo-EM structures and proof of concept data for cryo-EM using your own samples.

- ▶ [Find facilities offering cryo-EM instrument access for COVID-19 related research](#)
- ▶ [Find facilities offering cryo-EM instrument access for academics](#)

Ready to bring the power of cryo-EM into your own lab?

Thermo Fisher Scientific offers the expertise to support both your investment and your research endeavors. We can help you [precisely plan](#) every aspect of your laboratory to control building costs and limit downtime. Our [cryo-EM application specialists](#) can assist with everything from sample preparation to data processing. We provide [long-term ongoing support](#), as well as [flexible financing options](#) to get you in the game.

THERMO FISHER SCIENTIFIC

Thermo Fisher Scientific is the world leader in serving science. Our mission is to enable our customers to make the world healthier, cleaner, and safer. We serve life science researchers from academic institutions and the pharmaceutical industry, providing an unmatched combination of complete workflow solutions ranging from cryo-EM structural determination of macromolecular complexes and protein sociology in the native state to reconstruction of 3D architecture of tissues and cells. Our solutions help researchers unlock the mysteries of underlying protein function and cellular processes and bridge the gap between basic scientific research to design of new therapies.

REFERENCES

All structures made with ePMV. G.T. Johnson et al., "ePMV embeds molecular modeling into professional animation software environments," *Structure*, 19:293-303, 2011.

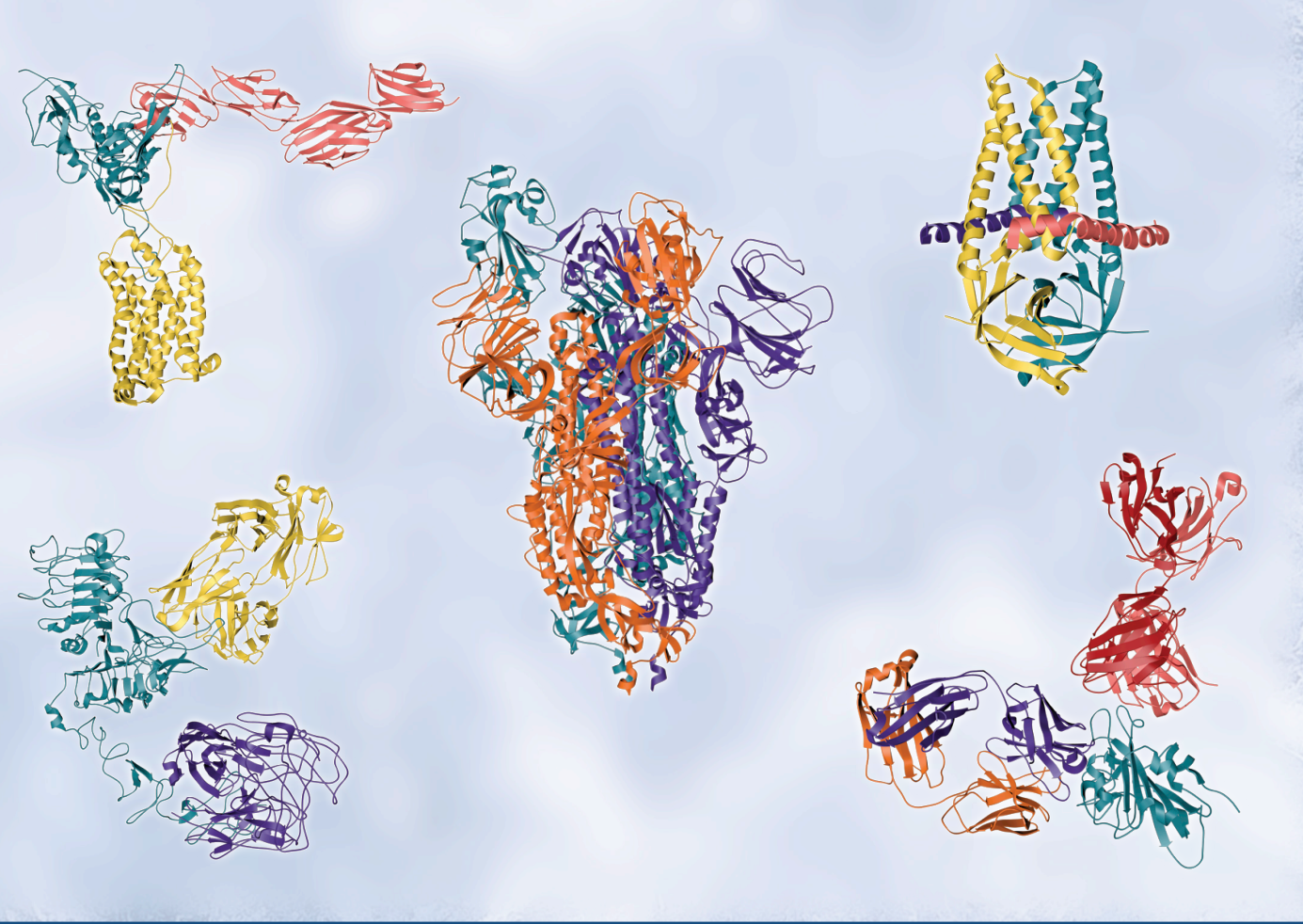
1. J. Dubochet, A.W. McDowell, "Vitrification of pure water for electron microscopy," *J Microsc*, 124:3-4, 1981.
2. D. Wrapp et al., "Cryo-EM structure of the 2019-nCoV spike in the prefusion conformation," *Science*, 367:1260-63, 2020.
3. J. Hansen et al., "Studies in humanized mice and convalescent humans yield a SARS-CoV-2 antibody cocktail," *Science*, 369:1010-14, 2020.
4. S. Ghosh et al., "β-coronaviruses use lysosomes for egress instead of the biosynthetic secretory pathway," *Cell*, 183:1520-35.e14, 2020.
5. C.-M. Chan et al., "The ion channel activity of the SARS-coronavirus 3a protein is linked to its pro-apoptotic function," *Int J Biochem Cell Biol*, 41:2232-39, 2009.
6. C. Castaño-Rodríguez et al., "Role of severe acute respiratory syndrome coronavirus viroporins E, 3a, and 8a in replication and pathogenesis," *mBio*, 9, 2018.
7. D.M. Kern et al., "Cryo-EM structure of the SARS-CoV-2 3a ion channel in lipid nanodiscs," *bioRxiv*, 2021.
8. G. Zhao et al., "Mature HIV-1 capsid structure by cryo-electron microscopy and all-atom molecular dynamics," *Nature*, 497:643-46, 2013.
9. M.M. Shaik et al., "Structural basis of coreceptor recognition by HIV-1 envelope spike," *Nature*, 565:318-23, 2019.
10. Z.T. Berendsen et al., "Visualization of the HIV-1 Env glycan shield across scales," *PNAS*, 117:28014-25, 2020.
11. C.K. Deniston et al., "Structure of LRRK2 in Parkinson's disease and model for microtubule interaction," *Nature*, 588:344-49, 2020.
12. R. Watanabe et al., "The in situ structure of Parkinson's disease-linked LRRK2," *Cell*, 182:1508-18.e16, 2020.
13. D. Nutt et al., "Psychiatric psychiatry's brave new world," *Cell*, 181:24-28, 2020.
14. K. Kim et al., "Structure of a hallucinogen-activated Gq-coupled 5-HT2A serotonin receptor," *Cell*, 182:1574-1588.e19, 2020.
15. M.D. Galbraith et al., "Therapeutic targeting of transcriptional cyclin-dependent kinases," *Transcription*, 10:118-36, 2019.
16. B.J. Greber et al., "The cryo-electron microscopy structure of the human CDK-activating kinase," *PNAS*, 117:22849-57, 2020.
17. N. Iqbal, N. Iqbal, "Human epidermal growth factor receptor 2 (HER2) in cancers: overexpression and therapeutic implications," *Mol Biol Int*, 2014:852748, 2014.
18. Y. Hao et al., "Cryo-EM structure of HER2-trastuzumab-pertuzumab complex," *PLOS ONE*, 14:e0216095, 2019.
19. E. Nogales, "The development of cryo-EM into a mainstream structural biology technique," *Nat Methods*, 13:24-27, 2016.
20. S.H.W. Scheres et al., "Modeling experimental image formation for likelihood-based classification of electron microscopy data," *Structure*, 15:1167-77, 2007.
21. K.M. Yip et al., "Breaking the next cryo-EM resolution barrier - atomic resolution determination of proteins!" *bioRxiv*, 2020.
22. T. Nakane et al., "Single-particle cryo-EM at atomic resolution," *bioRxiv*, 2020.

TheScientist
EXPLORING LIFE, INSPIRING INNOVATION

ThermoFisher
SCIENTIFIC

SOLVING STRUCTURES IN A FLASH

SCIENTIFIC BREAKTHROUGHS WITH CRYOGENIC ELECTRON MICROSCOPY



TheScientist
EXPLORING LIFE, INSPIRING INNOVATION

ThermoFisher
SCIENTIFIC

Tundra Cryo-TEM

Cryo-TEM that is cost effective and easy to use, bringing cryo-electron microscopy to every biochemistry laboratory.



The Thermo Scientific™ Tundra™ Cryo-Transmission Electron Microscope (Cryo-TEM) is a dedicated structure-determination solution designed to bring single particle analysis to every biochemistry laboratory. The Tundra Cryo-TEM enables fresh possibilities in structural research and analysis that have previously been out of reach thanks to simplified sample loading and built-in AI that makes it easier for beginners to use while also increasing productivity for experts.

- It is easier to use than typical cryo-TEM instruments
- Fits into a standard lab space
- Matches grant mechanisms and funding opportunities globally

The Tundra Cryo-TEM is a powerful tool that can help answer your most challenging research questions, offering structural determination at biologically relevant resolutions.

▶ [Find out more at thermofisher.com/Tundra](#)

