

Fragment-Based Drug Discovery

The long and difficult journey of drug discovery depends greatly on the quality of lead compounds. High Throughput Screening (HTS) has been heavily relied upon to search for such lead compounds. Recently however, Fragment-Based Drug Discovery (FBDD), which is a relatively new technology to identify potential fragments under 300 Da and optimize them into drugs, has been a topic of much focus in the field. In 1996, Fesik introduced FBDD as an alternative approach to HTS for the first time in the literature¹.

One desirable quality of a compound library is high completeness, in other words large chemical space occupancy². Chemical space is a concept in cheminformatics referring to the three-dimensional space constructed with predicted virtual molecules made up of the chemical elements represented, their number, and combination. The smaller the compounds that make up the library, the more comprehensive the chemical library can be. This means that the probability of finding effective compounds is higher in a fragment library consisting of small compounds.

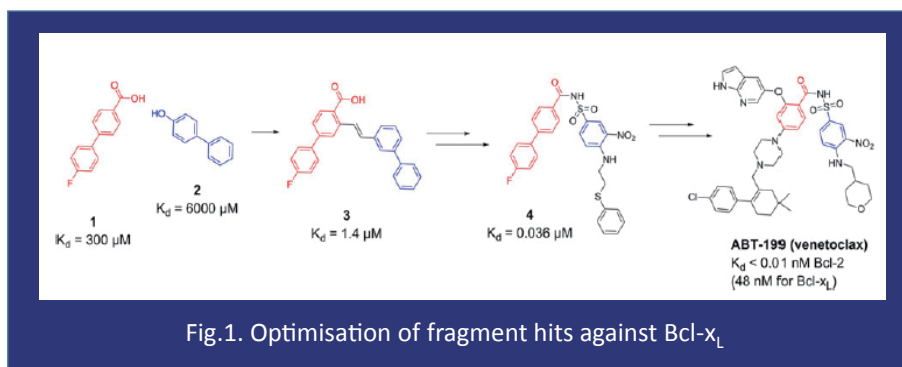
Hann reported that small and simple compounds have a lower probability of non-specific interaction with proteins and a higher probability of specific interaction with a target protein, compared to large and complex compounds³. This implies a high chance of success using fragment screening. Small lead fragments generally fulfill the criteria of Lipinski's "Rule of Five", thus lending themselves to the ability to expand lead fragments synthetically with a high degree of freedom while maintaining desirable drug-like physical and chemical properties.

However, the physiological inhibitory activity of small fragments are too weak to detect using conventional assay methods like EL

ISA. To identify binding fragments, it is often necessary to combine a number of biophysical measurement systems such as NMR, X-ray crystallography, thermal shift assay, and SPR, thus obtaining diversified information. The throughput of these techniques is inherently low.

Venetoclax, a Bcl-xL inhibitor launched as a treatment for chronic lymphocytic leukemia, is a successful example of a drug discovery program that started from fragments. Venetoclax was created by connecting two fragments (Fragment 1 and Fragment 2 in Fig. 1) discovered by FBDD with a linker, then optimizing them⁴. The K_d values of Fragment 1 and Fragment 2 were several hundred to several thousand μM , but the K_d value of Venetoclax was improved to 48 nM.

Carna is excited to introduce the benefits of unique and proprietary Weak Affinity Chromatography (WACTM) fragment screening services to its customers. This methodology is based on the combination of passing libraries of fragments over a protein covalently immobilized on a standard high-performance liquid chromatography (HPLC) column, and mass spectrometry. The K_D value is directly calculated from the difference of retention time between the column with immobilized protein and the control, empty column (Fig. 2). Since the fragments eluted from the column are directly detected, the throughput is high and it is possible to measure 3,000-4,000 fragments per week. The results



of WAC are reported to highly correlate with those from NMR and SPR, at 88% and 83%, respectively⁵. WAC makes it possible to identify fragments with a wide range of affinity, from 0.1 μ M to 10mM, and decreases the necessary sample protein amount to 1/5-1/10 of other methods, while at the same time having a lower cost than SPR or X-ray crystallography. Services are offered using our 1,300-fragment library for full screening, or you can have the flexibility to utilize your own fragment library.

Please don't hesitate to [contact us](#) for further information or if you have questions.

• **Weak Affinity Chromatography (WAC™) Fragment Screening**

References :

- 1) Science. 1996; 274(5292):1531-4. Fesik SW.
- 2) Yakugaku Zasshi, 2010; 130(3) 315-323. Tanaka D.
- 3) J Chem Inf Comput Sci. 2001; 41(3):856-64. Hann MM.
- 4) Essays Biochem. 2017; 61(5):475-484. Price AJ.
- 5) Anal Chem. 2013; 85(14):6756-66. Meiby E.

