
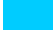



# 2025 Schedule for Carna's Profiling Services

The new service requests will be launched every Monday.

Please ensure that your complete set of test samples and the application form are received by us by **16:00 JST of each cutoff date** highlighted in yellow below.

For your Preincubation study request, please contact us at [info@carnabio.com](mailto:info@carnabio.com)

	: Test sample cutoff dates for next assay performance
	: %Inhibition Assays in MSA / IMAP <sup>®</sup> / ADP-Glo <sup>™</sup>
	: IC50 Assays in MSA / IMAP <sup>®</sup> / ADP-Glo <sup>™</sup>
Date in Red	: Off days

June						
Sun	Mon	Tue	Wed	Thu	Fri	Sat
1	2	3	4	5	6	7
8	9	10	11	12	13	14
15	16	17	18	19	20	21
22	23	24	25	26	27	28
29	30					

July						
Sun	Mon	Tue	Wed	Thu	Fri	Sat
		1	2	3	4	5
6	7	8	9	10	11	12
13	14	15	16	17	18	19
20	21	22	23	24	25	26
27	28	29	30	31		

August						
Sun	Mon	Tue	Wed	Thu	Fri	Sat
					1	2
3	4	5	6	7	8	9
10	11	12	13	14	15	16
17	18	19	20	21	22	23
24	25	26	27	28	29	30
31						

September						
Sun	Mon	Tue	Wed	Thu	Fri	Sat
	1	2	3	4	5	6
7	8	9	10	11	12	13
14	15	16	17	18	19	20
21	22	23	24	25	26	27
28	29	30				

October						
Sun	Mon	Tue	Wed	Thu	Fri	Sat
			1	2	3	4
5	6	7	8	9	10	11
12	13	14	15	16	17	18
19	20	21	22	23	24	25
26	27	28	29	30	31	

November						
Sun	Mon	Tue	Wed	Thu	Fri	Sat
						1
2	3	4	5	6	7	8
9	10	11	12	13	14	15
16	17	18	19	20	21	22
23	24	25	26	27	28	29
30						

December						
Sun	Mon	Tue	Wed	Thu	Fri	Sat
	1	2	3	4	5	6
7	8	9	10	11	12	13
14	15	16	17	18	19	20
21	22	23	24	25	26	27
28	29	30	31			

Question? Please feel free to contact us:

As of April 1, 2025

Carna Biosciences, Inc.  
 BMA 3F, 1-5-5 Minatojima-Minami, Chuo, Kobe 650-0047 JAPAN  
 E-mail: [info@carnabio.com](mailto:info@carnabio.com)  
 TEL: +81 78-302-7091

[www.carnabio.com](http://www.carnabio.com)

Note: Please visit our website for the latest schedule. Information may be changed without notice.

© 2012 Carna Biosciences, Inc.