

<b>Company</b>	
<b>Name</b>	
<b>Address</b>	
<b>Phone</b>	
<b>e-mail</b>	

**Short description of project, relevance to Covid 19**

--

		Test Item		
		1	2	3
<b>Therapeutic agent</b>	<b>Small molecule</b>			
	<b>Oligonucleotide</b>			
	<i>RNA</i>			
	<i>siRNA</i>			
	<i>mRNA</i>			
	<i>other (please describe)</i>			
	<b>Antibody</b>			
	<b>Vaccine</b>			
	<i>Protein</i>			
	<i>Peptide</i>			
	<i>Oligosaccharide</i>			
	<b>Cytokine</b>			
	<b>Other (please describe)</b>			

<b>Preparations</b>	Solvent			
	Concentration (=200x max. dosing concentration)			
	Number of vials			
	Volume			
	Suggested serial dilution			
	Storage recommendations			

<b>Expected compound delivery date</b>	
--	--

<b>Signatures &amp; stamp</b>			
	Principle Investigator	Department Manager	Company stamp

**Please fill out, sign and e-mail to [info@insphero.com](mailto:info@insphero.com) to fight Covid-19 together!**