

FFPE BLOCK DATA



Item Number	Batch Number	Sample Type	Number Samples Available	HPL	Barcode	Sample Harvest Date (Collection Date)	Species	Donor Sex	Donor Age	Donor Race	Tissue	Disease State
45048	818019	FFPE block	1	15000	HUZ652451 0400	1/14/2012	Human	F	18	More than one race	Placenta	Normal
45048	818020	FFPE block	1	15001	HUZ652440 0400	1/15/2012	Human	F	18	Black or African American	Placenta	Normal
45048	818021	FFPE block	1	15177	HUZ659843 0400	5/4/2012	Human	F	30	More than one race	Placenta	Normal
45048	818022	FFPE block	1	15297	HUZ659703 0400	8/1/2012	Human	F	20	Black or African American	Placenta	Normal
45048	818023	FFPE block	1	15306	HUZ659484 0401	8/8/2012	Human	F	19	Black or African American	Placenta	Normal
45233	818024	FFPE block	1	14982	HUZ652341 0408	1/2/2012	Human	F	21	Black or African American	Placenta	Preeclampsia
45233	818025	FFPE block	1	14988	HUZ652473 0408	1/6/2012	Human	F	20	Black or African American	Placenta	Preeclampsia
45233	818026	FFPE block	1	15019	HUZ657961 0408	1/22/2012	Human	F	21	Unknown	Placenta	Preeclampsia
45233	818028	FFPE block	1	14984	HUZ652086 0408	1/4/2012	Human	F	24	Black or African American	Placenta	Preeclampsia
45233	818030	FFPE block	1	15009	HUZ652263 0408	1/18/2012	Human	F	27	Black or African American	Placenta	Preeclampsia
45234	818031	FFPE block	1	14990	HUZ652325 0408	1/9/2012	Human	F	35	Black or African American	Placenta	Gestational Diabetes
45234	818032	FFPE block	1	15103	HUZ658036 0408	3/12/2012	Human	F	37	Black or African American	Placenta	Gestational Diabetes
45234	818033	FFPE block	1	15064	HUZ658012 0408	2/18/2012	Human	F	22	Black or African American	Placenta	Gestational Diabetes
45234	818034	FFPE block	1	15105	HUZ658201 0408	3/14/2012	Human	F	24	White	Placenta	Gestational Diabetes
45234	818035	FFPE block	1	15165	HUZ659790 0405	4/23/2012	Human	F	27	Black or African American	Placenta	Gestational Diabetes
45235	818036	FFPE block	1	16059	HUZ667836 0407	6/28/2013	Human	F	28	Black or African American	Placenta	Basal Plate, Normal
45235	818037	FFPE block	1	16159	HUZ668041 0406	6/24/2013	Human	F	25	Black or African American	Placenta	Basal Plate, Normal
45235	818038	FFPE block	1	16054	HUZ667894 0407	6/25/2013	Human	F	29	Black or African American	Placenta	Basal Plate, Normal
45235	818040	FFPE block	1	16114	HUZ667834 0406	7/19/2013	Human	F	32	Black or African American	Placenta	Basal Plate, Normal
45235	818042	FFPE block	1	16133	HUZ667756 0406	7/25/2013	Human	F	21	Black or African American	Placenta	Basal Plate, Normal