

Home Genomes Blat Tables Gene Sorter PCR Session FAQ Help

### Human (*Homo sapiens*) Genome Browser Gateway

The UCSC Genome Browser was created by the [Genome Bioinformatics Group of UC Santa Cruz](#).  
Software Copyright (c) The Regents of the University of California. All rights reserved.

clade genome assembly position or search term image width  
Mammal Human Mar. 2006 chr2:1-242,951,149 800 submit

[Click here to reset](#) the browser user interface settings to their defaults.

add custom tracks  clear position

Add custom tracks をクリック

#### About the Human Mar. 2006 (hg18) assembly (sequences)

The March 2006 human reference sequence (NCBI Build 36.1) was produced by the International Human Genome Sequencing Consortium.

#### Sample position queries

A genome position can be specified by the accession number of a sequenced genomic clone, an cDNA, an EST, an STS, a maplet, or a cytological band, a chromosomal coordinate, a maplet, or a cytological band.



Home Genomes Genome Browser Blat Tables Gene Sorter PCR Session FAQ Help

### Add Custom Tracks

clade Mammal genome Human assembly Mar. 2006 [hg18]

Display your own data as custom annotation tracks in the browser. Data must be formatted in [BED](#), [bigBed](#), [BEDGRAPH](#), [GFF](#), [GTF](#), [WIG](#), [bigWig](#), [MAF](#) or [PSL](#) formats. To configure the display, set [track](#) and [browser](#) line attributes as described in the [User's Guide](#). URLs for data in the bigBed and bigWig formats must be embedded in a track line in the box below. Publicly available custom tracks are listed [here](#). Examples are [here](#).

Paste URLs or data: Or upload:

Optional track documentation: Or upload:

参照をクリックし、bedファイルを選択後、Submitをクリック

Home Genomes Genome Browser Blat Tables Gene Sorter

### Manage Custom Tracks

genome: Human assembly: Mar. 2006 [hg18]

Name	Description	Type	Doc	Items	Pos	delete
<a href="#">User Track</a>	User Supplied Track	bed		1109682	chr2:	<input type="checkbox"/>

add custom tracks  
 go to genome browser  
 go to table browser

先ほど選択したbedファイルが trackとして組み込まれました

Go to genome browserをクリック



Home Genomes Blat Tables Gene Sorter PCR DNA Convert Ensembl NCBI PDF/PS Session Help

## UCSC Genome Browser on Human Mar. 2006 Assembly (hg18)

move <<< << < > >> >>> zoom in 1.5x 3x 10x base zoom out 1.5x 3x 10x

position/search chr2:20,000,000-20,100,000 jump clear size 100,001 bp. configure

chr2 (p24.1) 21 14 p12 3435

Scale 50 kb  
 chr2: 20010000| 20020000| 20030000| 20040000| 20050000| 20060000| 20070000| 20080000| 20090000|

User Track  
 User Supplied Track  
 Chromosome Bands Localized by FISH Mapping Clones

Human mRNAs from GenBank  
 Human ESTs That Have Been Spliced  
 Agilent Microarray Probesets

Ag CGH 1x1m  
 Ag HRD 1x1m  
 Ag CGH 2x400k  
 Ag CGH 4x180k  
 Ag CGH 8x60k  
 Ag CNV 2x400k  
 Ag CGH 1x244k  
 Ag CGH 2x105k  
 Ag CNV 2x105k  
 Ag CGH 4x44k

Vertebrate Multiz Alignment & Conservation (44 Species)  
 Placental Mammal Basewise Conservation by PhyloP

Mammal Cons

User Track: 黒いところがプローブの位置を示しています。

Affy SNP 6.0  
Affy SNP 6.0 SV  
Illumina IM-Duo  
RepeatMasker

move start Click on a feature for details. Click or drag in the base bars on left for track options and descriptions.

< 2.0 >

default tracks hide all manage custom tr

collapse all expand all

Use drop-down controls below and press refresh to alter tracks displayed.  
Tracks with lots of items will automatically be displayed in more compact modes.

User Trackの項目を"Full"にすると、位置が重なっているものが別々に表示されます。

Home Genomes Blat Tables Gene Sorter PCR DNA Convert Ensembl NCBI PDF/PS Session Help

UCSC Genome Browser on Human Mar. 2006 Assembly (hg18)

move <<< << < > >> >>> zoom in 1.5x 3x 10x base zoom out 1.5x 3x 10x

position/search chr2:20,000,000-20,100,000 jump clear size 100,001 bp. configure

chr2 (p24.1) 21 14 12 34 35

Scale chr2: 50 kb  
20010000 20020000 20030000 20040000 20050000 20060000 20070000 20080000 20090000

User Supplied Track

Human mRNAs From GenBank  
Spliced ESTs  
Human ESTs That Have Been Spliced  
Agilent Microarray Probesets

Ag CGH 1x1M  
Ag HRD 1x1M  
Ag CGH 2x400K  
Ag CGH 4x180K  
Ag CGH 3x50K

Home Genomes Blat Tables Gene Sorter PCR DNA Convert Ensembl NCBI PDF/PS Session Help

UCSC Genome Browser on Human Mar. 2006 Assembly (hg18)

move <<< << < > >> >>> zoom in 1.5x 3x 10x base zoom out 1.5x 3x 10x

position/search chr2:20,000,000-20,001,000 jump clear size 1,001 bp. configure

chr2 (p24.1) 21 14 12 34 35

Scale chr2: 500 bases  
20000100 20000200 20000300 20000400 20000500 20000600 20000700 20000800 20000900

User Supplied Track

Human mRNAs From GenBank

Chromosome Bands Localized by FISH Mapping Clones

A\_16\_F57819884  
A\_16\_F15575485  
A\_16\_F09327341  
A\_18\_F21492694  
A\_16\_F15575485  
A\_16\_F09327342  
A\_18\_F21491820  
A\_16\_F15575487  
A\_18\_F21494492  
A\_16\_F21492216  
A\_16\_F15575488  
A\_18\_F21492727  
A\_18\_F13139733  
A\_16\_F30541630  
A\_16\_F30541631  
A\_16\_F30541632  
A\_16\_F30541633  
A\_16\_F30541634  
A\_16\_F30541635

さらにzoom-inすると~60merのプローブの位置がより明確に分かります。